

Part III, Endotoxin Test Concerns of Biologics: *LER From a Broad Biologics Test Perspective*

Kevin Williams

Senior Scientist, R&D-Endotoxin, Lonza Walkersville, Inc.

Overview

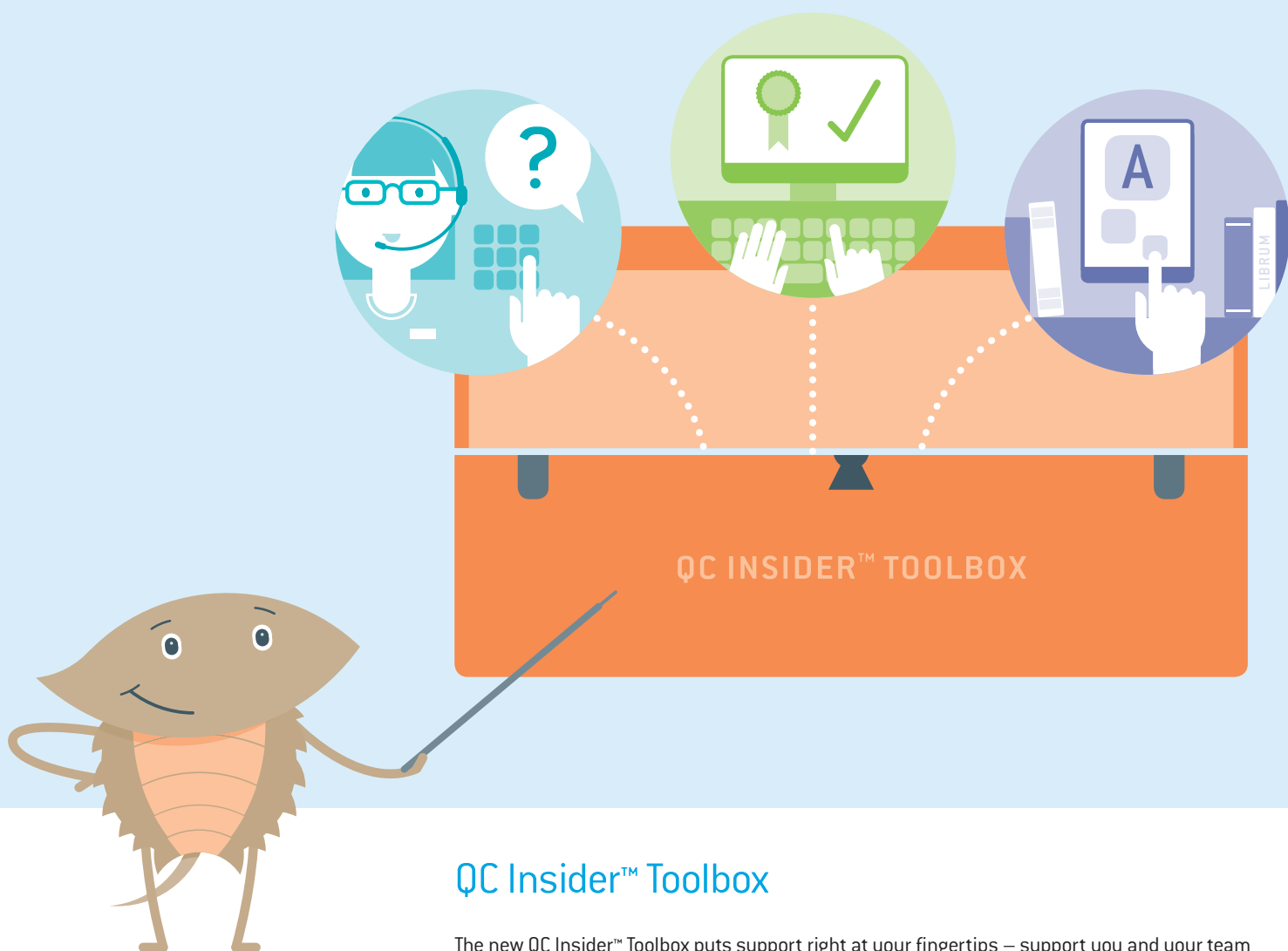
Recent conferences have demonstrated the polarizing nature of the Low Endotoxin Recovery debate, including those in Berlin (PDA Europe), Iselin, NJ (PMF Bacterial Endotoxin Summit), and Bethesda, MD (PDA Global Micro). The debate has centered on the significance of samples subjected to LER conditions, which includes polysorbate and a chelator (citrate or phosphate buffer). Samples subjected to these conditions undergo a masking effect that renders undiluted spike into hold-time studies impossible to recover without special treatment. Industry participants are debating the potential for harm from such samples, in the unlikely event they were contaminated by biologics manufacturing processes. Three points outline a broad biologics-test-based endotoxin contamination control perspective:

1. Historical endotoxin detoxification studies
2. The emerging view of endotoxin as an IIRMI
3. A perspective on merging of the current pyrogen and emerging IIRMI view

Historical Endotoxin Detoxification Studies

There is a large volume of knowledge derived from the search for adjuvant candidates to aid the efficacy of vaccines that dates back over 80 years. These searches focused on the idea that one can treat purified lipopolysaccharide or natural endotoxin with harsh, or not so harsh conditions (see Figure 1) and produce a changed LPS structure that is no longer pyrogenic. However, while not pyrogenic, the endotoxin subjected to detoxifying conditions maintains its ability to stimulate the immune system. This immune stimulation is exactly what vaccine manufacturers want in order to direct the body's immune system against clean, recombinant proteins that do not, of

Endotoxin Expertise At Your Fingertips™



SUPPORT
TRAINING
LIBRARY

QC Insider™ Toolbox

The new QC Insider™ Toolbox puts support right at your fingertips – support you and your team can access any time and from anywhere in the world.

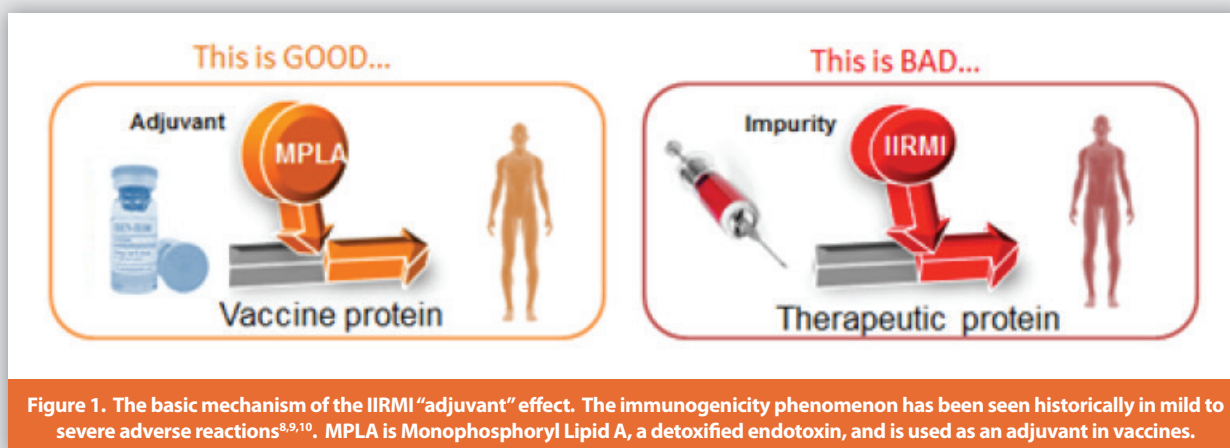
QC Insider™ Support – One-on-one guidance, software support, and more from our QC experts

QC Insider™ Training – Self-directed training resources that will help you build on your QC expertise

QC Insider™ Library – Product instructions and technical guides your team needs to succeed

Turn your endotoxin testing team into QC Insider™ experts
today by visiting www.lonza.com/qcinsider

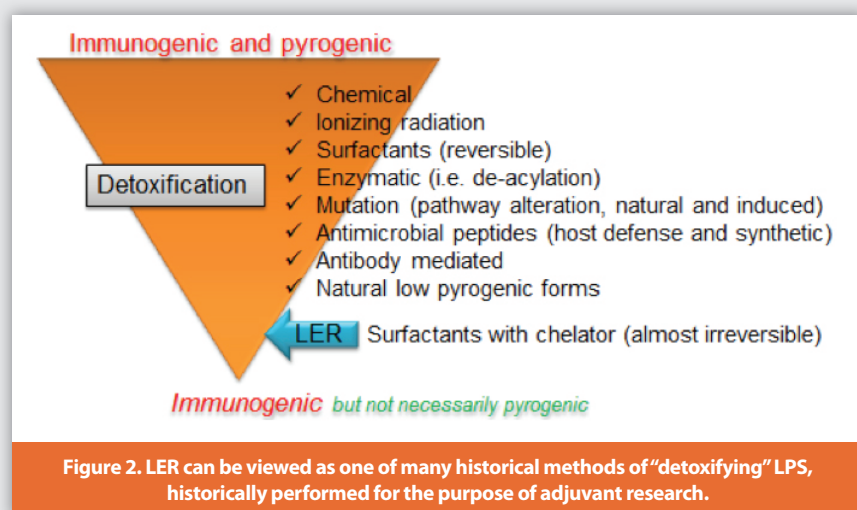
www.lonza.com



themselves, elicit an immune response¹. In this case, the stimulation of anti-drug antibodies (ADA) against recombinant drug proteins from bacterial and other antigens including Hepatitis B², Human papilloma virus³, various cancer vaccines^{4,5,6}, and Malaria⁷ is beneficial to the vaccine effort. The antigen-adjuvant effect forms the basis of vaccination, which is the most successful form of medical intervention ever employed. The downside of the so-called “adjuvant effect” involving clean proteins is that the same mechanism can serve to bring about ADAs toward life-saving therapeutic proteins via the inclusion of low levels of inadvertent microbial contaminants. While the LAL reduction associated with O-deacylation and dephosphorylation has

long been known, what is less recognized is that “detoxification” results in greatly muted LAL and rabbit pyrogen test activity (“when monophosphoryl lipid A’s and lipid X were similarly tested, they showed very low pyrogenicity”¹¹), however the adjuvant activity remains¹². A wide variety of historically accumulated methods of detoxification are shown in Figure 2. These include: chemical,^{13,14,15,16} ionizing radiation,^{17,18} use of surfactants¹⁹ (reversible), enzymatic^{20,21} mutation^{22,23,24} (natural and induced), antimicrobial peptides²⁵, antibody mediated²⁵ natural low pyrogenic forms²⁶ and LER. A review of practices that do not incinerate or completely remove the functional LPS PAMP might be in order from the “endotoxin as

IIRMI” view for biologics processing. Detoxification does not remove the adjuvant effect of MPLA but rather significantly diminishes the pro-inflammatory effect²⁷. This has been seen in methods used for depyrogenation as well, including Gamma irradiation, as seen in the treatment of *Salmonella* resulting in removal of its pyrogenicity, while allowing it to retain its immunogenicity-inducing capability²⁸. Similarly, irradiated LPS “retains the adjuvant activity of LPS, and it serves as a good adjuvant for inactivated virus vaccines.” LER solutions that are purposefully spiked or inadvertently contaminated could be considered a type of “detoxified” endotoxin. The search for compounds that utilize the immune stimulation properties of LPS without pro-inflammatory effects that include fever is ongoing, as many subunit vaccines do not have the ability to stimulate the immune system. In the realm of endotoxin testing, if one is singularly worried about the pyrogenicity of a sample, then it could come to play out that LER-subjected drug formulations are not particularly pyrogenic, although there is conflicting rabbit pyrogen data^{29,30}. However, if one is worried that a given LER-prone protein formulation could increase therapeutic protein immunogenicity if such LPS monomers are present, then one would want to detect and preclude the presence of endotoxin monomers or otherwise “detoxified” endotoxin solutions that retain the potential to be recognized by mammalian immune systems.



Achieving new standards in performance takes
years of innovative R&D. This is the result.



Vcaps® Plus hypromellose capsules
for pharmaceutical products – true pH and ionic media independence
in disintegration for optimized delivery and performance.

Some HPMC capsules are formulated with secondary gelling agents and ionic gel promoters – which may slow down release of active ingredients. Capsugel's Vcaps® Plus HPMC capsules are made through a proprietary manufacturing process that eliminates the need for gelling systems. They disintegrate and release their contents independent of pH and ionic strength of the test media. To discover how Vcaps® Plus capsules can optimize performance, improve product stability, and reduce both variability and timelines in your drug development, download our free whitepaper.

VcapsPlus.com 888-783-6361

CAPSUGEL®

© 2015 Capsugel Belgium NV All rights reserved.

“

There is clearly concern here centered on a level of bacterial endotoxin that is below the level associated with the ability to bring about fever.

”

An Emerging View of Endotoxin as an IIRMI

An emerging view of endotoxin as a contaminant comes from FDA laboratory studies (Verthelyi and Wang³¹; Haile et al.³²) that demonstrated that low levels of contaminants including LPS that they have termed IIRMIs or “innate immune response modulating impurities” can, either alone or synergistically to greater effect, stimulate the immune system against therapeutic proteins at a level that may be well below the level considered necessary for pyrogen or endotoxin preclusion testing (5 EU/kg). Studies in mice showed that LPS and bacterial DNA added to a recombinant protein (r-Erythropoietin) served to “break tolerance to self”.

Researchers have shown both *in vitro* and *in vivo* that synergistically IIRMI's are active at lower levels than when present alone³¹:

This synergistic effect was then confirmed *in vivo*, as studies showed that the combination of 10 ng of LPS and 500 ng of CpG ODN, which do not induce an immune response when present individually, were sufficient to promote the immunogenicity of proteins and contribute to a clinically relevant break in tolerance to self.

They showed that mice given r- Erythropoietin had mild transient side effects, while mice subject to r- Erythropoietin with low levels of LPS and bacterial DNA contaminant developed long-lasting anemia. Thus, the attack on the r- Erythropoietin turned into a full-fledged attack on the mice's own natural proteins. This is among the worst forms of undesirable immune reactions. Furthermore, either of the components acting alone (LPS or DNA) at very low levels did not show an adverse effect. The levels involved for mice do not translate directly to humans, as mice are notoriously less affected by endotoxin relative to humans³³.

The IIRMI view is referred to as “emerging” here, but it is already included in the 2014 FDA Guidance Document: *Immunogenicity Assessment for Therapeutic Protein Products* in section 5, page 18, “Impurities with Adjuvant Activity”. While considered “non-binding,” the non-binding designation also applies to the current Q&A Guidance document. An excerpt from the former is shown below:

Adjuvant activity can arise through multiple mechanisms, including the presence of microbial or host-cell-related impurities

in therapeutic protein products (Verthelyi and Wang 2010; Rhee et al. 2011; Eon-Duval et al. 2012; Kwissa et al. 2012). These innate immune response modulating impurities (IIRMIs), including lipopolysaccharide, β -glucan and flagellin, high-mobility group protein B1 (HMGB1), and nucleic acids exert immune-enhancing activity by binding to and signaling through toll-like receptors or other pattern-recognition receptors present on B-cells, dendritic cells, and other antigen-presenting cell populations (Iwasaki and Medzhitov 2010; Verthelyi and Wang 2010). This signaling prompts maturation of antigen-presenting cells and/or serves to directly stimulate B-cell antibody production.

The researchers quoted by the Immunogenicity Guidance document (Verthelyi and Wang) describe the current conundrum in endotoxin detection. They discuss the relevant levels of endotoxin viewed as an IIRMI to those standardized for testing of pyrogens, by either *Limulus*-based methods (LAL and rFC) or rabbit pyrogen tests (RPT).

Of note, the current guidelines for setting limits on these impurities are not based upon their potential impact on product immunogenicity. For example, the current recommendations for endotoxin content in parenteral products (0.5 EU/kg/hr)³⁴ is based on its pyrogenic potential, while WHO recommendations for DNA content (<10 ng DNA/dose) are based on minimizing the risk of DNA integration (USP <85> and [45])³⁵...the levels of agonists sufficient to stimulate an innate immune response can be lower when multiple receptors are engaged.

In terms of immunogenicity³⁶, biologics have become increasingly safe over time with the realization that natural animal-derived proteins may be recognized as non-self, and improved manufacturing methods for recombinant proteins has lessened potential protein aggregation to further limit immunogenic reactions. The “humanization” of previously animal-based and chimeric monoclonal antibodies has also lowered immunogenicity rates. But a low level (and sometimes not so low level) of persistent proclivity toward immunogenicity remains and can be seen in clinical studies and marketed package inserts.

A Perspective on Merging of the Current Pyrogen and Emerging IIRMI View

There is a “trend” that could be interpreted as a rudimentary, perhaps utilitarian effort to merge the pyrogen and IIRMI preclusion concepts. How else is one to interpret the continued reduction of endotoxin tolerance limits (TL) associated with biologics? This can be seen in (a) Biologic License Application (BLA) requests to lower limits calculated by the TL = K/M method, (b) hold-time study recoveries from undiluted drug process and final solutions, and (c) the Q&A verbiage that makes it clear that current expectations are that testing be performed as low as possible after interference has been overcome. There is clearly concern here centered on a level of bacterial endotoxin that is below the level associated with the ability to bring about fever. Historically, one would calculate the tolerance limit (TL) by dividing the threshold pyrogenic response constant (K=5 EU/kg/hr) by the drug dose in

active dosage units per kilograms of body weight. The dose of a given biologic drug appears to be less and less a consideration since the advent of biologics. Today, the question becomes “How far below the TL should be validated and routinely tested?”

Alternatively, the concerns could also stem from the large number of additional solutions going in with biologics such as pre-medications for expected fever reactions (steroids, antipyretic, antihistamine, etc.), which are themselves (the biologic and pre-medications) typically each contained within large volume parenteral solutions (LVP). Note that the limit for LVPs is less stringent than other drug types (0.5 EU/mL). According to the IIRMI or adjuvant view, a lesser quality standard for a drug to be co-administered or subsequently administered with a biologics could contribute IIRMI. One wonders if a quality designation of “For use in Biologics” might serve to safeguard against emerging concerns associated with the adjuvant effect that can come from non-biologic sources when administered with biologics. This concern extends to potential contaminants added by compound pharmacy activities³⁷, which have recently come under fire and to which the FDA has responded with a new draft guideline³⁸.

Historically, there has been a singular focus on precluding the bacteria that produce pro-inflammatory, “endotoxic” endotoxins (*Enterobacteriaceae*). The revelation that non-pyrogenic endotoxin can be recognized for other effects³⁹ fits an underlying, longstanding theme that microbial artifacts injected into the blood stream may have significant effects that do not necessarily correlate with our ability to “see” them, analytically speaking, or correlate with their ability to produce fever. The mammalian physiological view of endotoxin is ultra-sophisticated when it comes to the detection of microbe invaders and their artifacts, given that it has accrued over a billion of years of evolution. The basic Lipid A PAMP should be viewed as a set of dials (phosphate, sugar, number and types of acyl chains-symmetrical/asymmetrical, substitutions, etc.) rather than an “on-off” button (pyrogenic or non-pyrogenic)⁴⁰. The activity of LPS at low levels is being borne out in studies of the low dose effect of endotoxin in various disease states (i.e. sepsis⁴¹, inflammation⁴², cancer⁴³, and cardiovascular disease⁴⁴).

In LER discussions, test users have wondered if endotoxins disassociated by LER conditions could pass along any harm to patients. The adjuvant model seems to answer this question as a detoxified, non-pyrogenic endotoxin such as FDA-approved adjuvant monophosphoryl lipid A (MPLA) is added to recombinant vaccine proteins to stimulate the immune system against vaccine proteins. While the detoxified form is not pyrogenic (detoxified), and users are testing for pyrogenic effects via rabbit pyrogen and bacterial endotoxin test methods instead of immunogenicity, it answers the question as to whether it is possible for such forms to potentially harm patients. It is not only possible, it is an effect routinely taken advantage of in the vaccine realm.

The difficulty LER raises for regulators and drug manufacturers is that the endotoxin test currently provides a multifunctional duty as a process control to exclude endotoxins. There is not another method or a better method to control endotoxins in processes other than by recognizing their presence by bacterial endotoxin testing. If the BET vantage is lost due to a common masking phenomenon (LER or other biologics’ inherent process conditions), then the presence

of pyrogen-causing and immunogenicity-causing endotoxins cannot be known. There have been developmental efforts to restore the activity lost by the LER phenomenon as shown by J. Chen of Genentech, J. Reich of Hyglos and here at Lonza. While demasking takes a little more time and labor and may not be perfect from the perspective of process control, it does allow a way forward that does not completely ignore the potential harm from endotoxin that would, if present as a masked contaminant, remain masked right up to the point of patient administration. Disaggregated endotoxin consists of the entire microbial PAMP and active endotoxic principle (lipid A) that has served for a billion years (plants) of metazoan recognition of prokaryotic invasion. There is now ample evidence that even non-pyrogenic endotoxin is immunogenic, and that this activity proceeds by a pathway separate from the pro-inflammatory pathway after the initial recognition via TLR4/MD-2 dimer complex^{45,46}.

Conclusion

The last thing biologics manufacturers intend is to include adjuvants in the use of therapeutic proteins by way of impurities. As we have seen, endotoxin adjuvants administered with vaccine proteins do indeed elicit effective, non-pyrogenic endotoxin responses. The need for an updated view is well stated by Haile, et al: “It is only the more recent understanding of the innate immune system’s biology that dictates the need of assessing a broader spectrum of known and unknown IIRMI in order to control or reduce the risk of unwanted immunogenicity by therapeutic proteins⁴⁷.” Modern biologics manufacturing concerns contrast with historic, purely pyrogen-centric activities that have represented an important but more minimal standard.

In conclusion, three posits may help form a working model towards a broader goal of containing endotoxin contamination by including the IIRMI perspective:

1. Endotoxin detection from a multipurpose perspective should be performed at levels lower than those dictated solely by the TL calculation.
2. Processes known to employ detoxifying conditions (such as LER) should pursue orthogonal test methods that supplement conventional methods.
3. The philosophical view that non-biologics cannot practically contribute to biologics contamination should be challenged given the adjuvant activity they can contribute to biologics via co-administration.

References

1. “A major complication of subunit vaccine development is that most recombinant proteins lack intrinsic immunostimulatory activity.” From: Putting endotoxin to work for us: monophosphoryl lipid A as a safe and effective vaccine adjuvant, C. R. Casella and T. C. Mitchell, Cell Mol Life Sci. 2008 October ; 65(20): 3231–3240.
2. Vaccine Adjuvant Systems containing monophosphoryl lipid A and QS-21 induce strong humoral and cellular immune responses against hepatitis B surface antigen which persist for at least 4 years after vaccination, G. Leroux-Roels et al., Vaccine 33 (2015) 1084–1091.

3. FDA Licensure of Bivalent Human Papillomavirus Vaccine (HPV2, Cervarix) for Use in Females and Updated HPV Vaccination Recommendations from the Advisory Committee on Immunization Practices (ACIP), Morbidity and Mortality Weekly Report (MMWR), May 28, 2010 / 59(20);626-629. Accessed on-line June 16, 2015.
4. Monophosphoryl Lipid A (MPL) as an Adjuvant for Anti-Cancer Vaccines: Clinical Results, Christopher W Cluff.
5. Christopher W. Cluff, Lipid A in Cancer Therapy, edited by Jean-François Jeannin. ©2009 Landes Bioscience and Springer Science+Business Media.
6. Effect of DETOX as an adjuvant for melanoma vaccine, Schultz, N. et al., Vaccine. 1995 Apr;13(5):503-8.
7. Randomized, double-blind, phase 2a trial of Falciparum Malaria vaccines RTS,S/AS01B and RTS,S/AS02A in malaria-naïve adults: safety, efficacy, and immunologic associates of protection, K.E. Kester et al., Jour. Infect. Dis., 2009; 200:337-46.
8. Drugs Highly Associated with Infusion Reactions Reported using Two Different Data-mining Methodologies, P.W. Moore, et al., J Blood Disorders Transf 2014, 5:2.
9. Cytokine Storm in a phase I trial of the anti-CD28 monoclonal antibody TGN1412, Suntharalingam, et al., N. Engl. J. Med., 355, pp. 1018-1028
10. Clinical review: Serious adverse events associated with the use of rituximab - a critical care Perspective, Kasi et al. Critical Care 2012, 16:231.
11. Influence of Fine Structure of Lipid A on Limulus Amebocyte Lysate Clotting and Toxic Activities, K. Takayama et al., Infection and Immunity, Aug. 1984, p. 350-355 Vol. 45, No. 2.
12. Development and Characterization of Synthetic Glucopyranosyl Lipid Adjuvant System as a Vaccine Adjuvant. Coler RN, Bertholet S, Moutafsi M, Guderian JA, Windish HP, et al. (2011) PLoS ONE 6(1): e16333.
13. The Immunobiology of TLR4 Agonists: From Endotoxin Tolerance to Immunoadjuvants, J. K. Bohannon et al., Shock. 2013 December; 40(6): 451-462.
14. Preparation and biological properties of a chemically modified Escherichia coli endotoxin of high immunogenic potency and low toxicity. J. Clin. Inv., 1961, 40, 1935.
15. The adjuvant monophosphoryl lipid A increases the function of antigen-presenting cells Int. De Becker, et al. Immunol, 2000.
16. Detoxification of Bacterial Endotoxin with Retention of Ability to Stimulate Non-Specific Resistance to Infection, H. Freedman et al., Exp Biol Med (Maywood) August 1961 107: 819-821.
17. Immunogenicity of irradiated Salmonella typhimurium cells and endotoxin, J.J. Preville, Jour. Bact. June 1968, 2165-2170.
18. Radio-detoxified endotoxin activates natural immunity: a review, F. József Jour. Pathophys. 2005.02.004.
19. Immunogenicity of Deoxycholate-disrupted Endotoxins, J. Bact. Jan. 1969, p. 13-15.
20. Crystal structure of LpxC, a zinc-dependent deacetylase essential for endotoxin biosynthesis, D. A. Whittington et al., Proc Natl Acad Sci U S A. 2003 Jul 8; 100(14): 8146-8150.
21. A Host Lipase Detoxifies Bacterial Lipopolysaccharides in the Liver and Spleen, B. Shao, et al., Jour. of Biol. Chem., VOL. 282, No. 18, May 4, 2007.
22. Bacterial outer membrane vesicles and vaccine applications, R. Acevedo et al., Front. Immunol., 24 March 2014.
23. Salmonella synthesizing 1-monophosphorylated LPS exhibits low endotoxic activity while retaining its immunogenicity, Jour. Immunol. 2011 Jul 1; 187(1): 412-423.
24. Detoxifying Escherichia coli for endotoxin-free production of recombinant proteins, U. Mamat et al., Microbial Cell Factories (2015) 14:57 (not immunogenic).
25. Parameters Involved in Antimicrobial and Endotoxin Detoxification Activities of Antimicrobial Peptides, Y. Rosenfeld et al., Biochemistry, 2008, 47 (24), pp 6468-6478.
26. Naturally occurring low biological reactivity lipopolysaccharides as vaccine adjuvants, Expert Rev. Vaccines, Darveau and Chilton, 12 (7), 707-709 (2013).
27. The Vaccine Adjuvant Monophosphoryl Lipid A as a TRIF-Biased Agonist of TLR4, V. Mata-Haro et al., Science 15, June 2007: Vol. 316 no. 5831 pp. 1628-1632.
28. Immunogenicity of irradiated Salmonella typhimurium cells and endotoxin, J.J. Preville, Jour. Bact. June 1968, 2165-2170.
29. Low endotoxin recovery: An FDA perspective, P.F. Hughes, et al., BioPharm Asia, published April 2015, accessed on-line June 12, 2015.
30. P. Hughes, "Biotech Manufacturing Assessment Branch, FDA/CDER," presentation at the PDA Conference (Bethesda, MD, Oct.21, 2014).
31. Trace Levels of Innate Immune Response Modulating Impurities (IIRMI) Synergize to Break Tolerance to Therapeutic Proteins. Verthelyi D., Wang V. Vij N, ed. PLoS ONE. 2010; 5(12):e15252.
32. Detection of Innate Immune Response Modulating Impurities in Therapeutic Proteins, L.A. Haile et al., PLOS ONE, DOI:10.1371/journal.pone.0125075, April 22, 2015. U.S. Department of Health and Human Services, Food and Drug Administration, Center for Drug Evaluation and Research (CDER), Center for Biologics Evaluation and Research (CBER), August 2014.
33. "Relative to the human response, mice are highly resilient to inflammatory challenge. For example, the lethal dose of endotoxin is 5-25 mg/kg for most strains of mice, whereas a dose that is 1,000,000-fold less (30 ng/kg) has been reported to cause shock in humans. Genomic responses in mouse models poorly mimic human inflammatory diseases." J. Seok, et al., Proc Natl Acad Sci U S A. 2013 Feb 26; 110(9):3507 -3512.
34. This value is 5 EU/kg/hr as the threshold pyrogenic response (K) via the parenteral route of administration.
35. Requirements for the use of animal cells as in vitro substances for the production of biologicals. 878. 8-9. 1998. World Health Organization. Requirements for Biological Substrates. No50.
36. In the context of Therapeutic Proteins product, immunogenicity refers to the immune response of the host against the Therapeutic Protein", The immunogenicity of therapeutic proteins- what you don't know can hurt YOU and the patient, J.A. Pedras-Vasconcelos, CMC and Immunogenicity Reviewer, Div. of Therapeutic Proteins, FDA, Fall 2014 (presentation accessed on-line June 22).
37. Compounding: Inspections, Recalls, and other Actions: <http://www.fda.gov/Drugs/GuidanceComplianceRegulatoryInformation/PharmacyCompounding/ucm339771.htm>
38. Mixing, Diluting, or Repackaging Biological Products Outside the Scope of an Approved Biologics License Application Guidance for Industry, DRAFT GUIDANCE, U.S. Department of Health and Human Services Food and Drug Administration Center for Drug Evaluation and Research (CDER) Center for Biologics Evaluation and Research (CBER) February 2015.
39. Insights into the species-specific TLR4 signaling mechanism in response to Rhodobacter sphaeroides lipid A detection, Anwar M.A. et al, Nature Scientific Reports, Dec. 2014, 5:7657.
40. Modulating the innate immune response by combinatorial engineering of endotoxin, Needham, B.D., et al, PNAS, Jan. 2013, Vol. 110, No. 4, 1464-1469.
41. Super-low Dose Endotoxin Pre-conditioning Exacerbates Sepsis Mortality, K. Chen et al., EBioMedicine 2 (2015) 324-333.
42. Molecular and cellular mechanisms responsible for cellular stress and low-grade inflammation induced by a super-low dose of endotoxin, B. Baker et al., J Biol Chem. 2014 Jun 6;289(23): 16262-16269
43. Therapeutic targeting of toll-like receptors for infectious and inflammatory diseases and cancer, O'Neill, L.A.J et al., Pharmacological Reviews, 2009, 61:177 -197
44. Low level bacterial endotoxin activates two distinct signaling pathways in human peripheral blood mononuclear cells, Blomkalns, A.L. et al., Journal of Inflammation 2011, 8:4.
45. Role of Adaptor TRIF in the MyD88-Independent Toll-Like Receptor Signaling Pathway, M. Yamamoto et al., Science, 1 Aug. 2003 Vol. 301.
46. Beyond MyD88 and TRIF pathways in Toll-like receptor signaling , V. Piras and K. Selvarajoo, Frontiers in Immunology, vol. 5, article 70, Feb. 2014.
47. Detection of Innate Immune Response Modulating Impurities in Therapeutic Proteins, L.A. Haile et al., PLOS ONE, DOI:10.1371/journal.pone.0125078, April 22, 2015.