

Controlling Variability to Increase Accuracy in the BET Assay

1

KAREN ZINK MCCULLOUGH
MMI ASSOCIATES

(908) 534-6463
karenzm@embarqmail.com

Definition of Accuracy

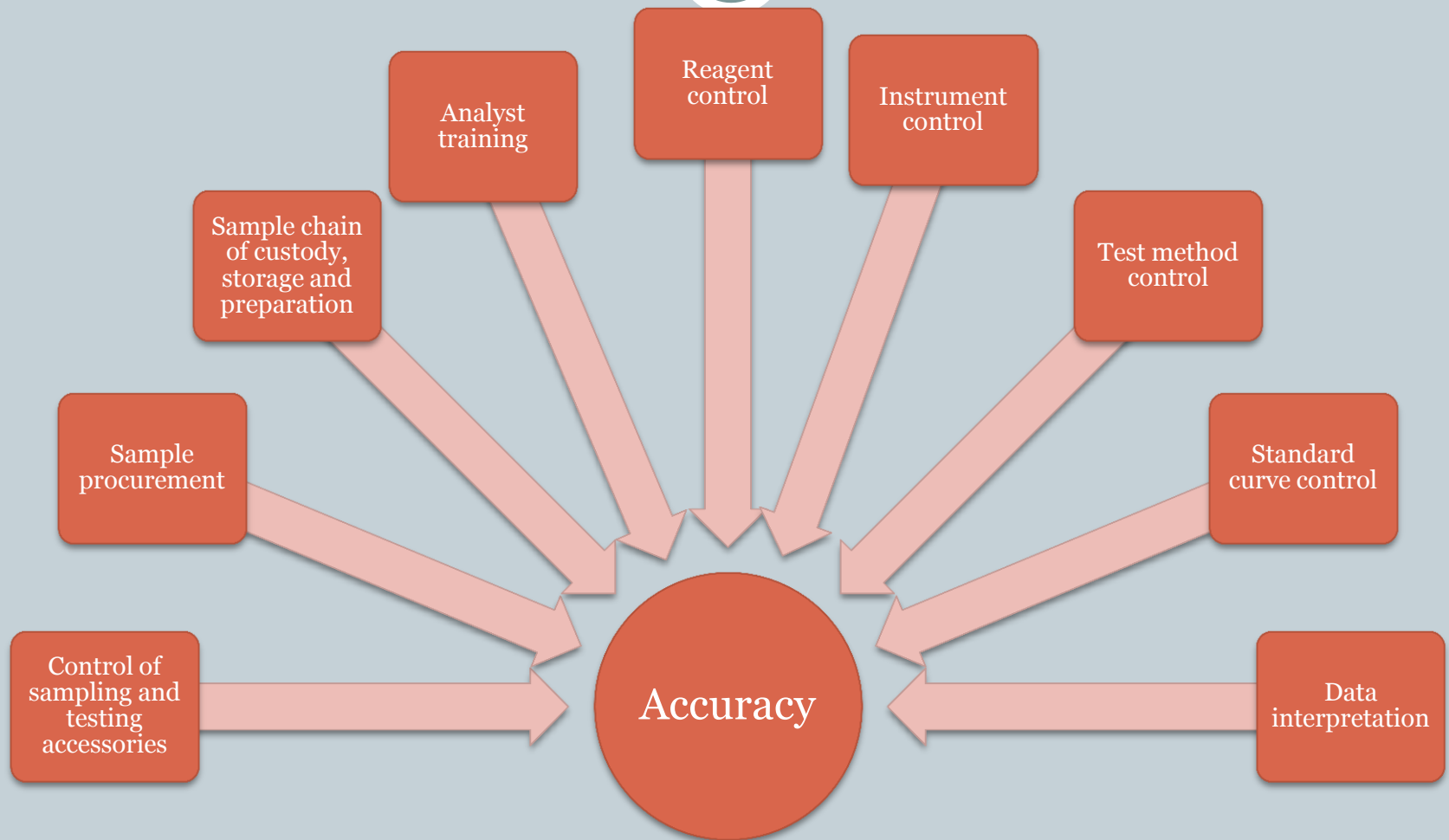
2

The degree to which the result of a measurement, calculation, or specification conforms to the correct value or a standard.

VARIABILITY in the BET assay affects accuracy

What Affects Accuracy in a BET Assay?

3



Control of Sampling and Testing Accessories

4

- Depyrogenated sampling equipment
 - Dry heat for glass bottles/metal scoops
 - Demonstration of lack of interference for plastics
- Testing accessories must be shown to be free of interference (plastics) – Reference USP <85>
 - Detectable endotoxin
 - Leachables
 - Assign a low limit and test according to AAMI ST72 or USP <161> (medical devices)
 - Polystyrene is usually OK. Polypropylene can pose problems
 - No requirement for spike/recovery in USP <85>, <161>, or AAMI ST72
 - Manufacturer's CoA should be confirmed

Sample Procurement

5

- Sample vessels are non-interfering
- Sample (water, in process product stream, raw materials) procured aseptically as not to introduce contamination
 - Clear training and supervision for samplers

Sample Chain of Custody, Storage, Preparation

6

- ❑ For non-steriles, transport in a manner that will not support microbial proliferation
 - Validate time/temperature of transport/hold time
- ❑ Validate storage conditions
- ❑ Always mix/vortex any sample prior to testing and between dilutions to assure that endotoxin is distributed
 - Especially important when working with CSE “spikes” because of the natural aggregation of the purified LPS
 - Important as well for natural endotoxin in samples, even though the surrounding cell wall components help to keep the endotoxin in solution

Analyst Training

7

- Assay is hugely analyst/technique dependent
- Need training in
 - Aseptic handling of samples during preparation
 - Preparing serial dilutions/special sample preparation
 - Pipetting
 - Serological and mechanical pipettors - technique
 - Analysis of data – when are tests invalid and when are they OOS?
- Best to have a combination of
 - “Book” training (theory and practical matters)
 - SOP training
 - Demonstration of competency

Reagent Control

8

- ❑ Confirm sensitivity of reagents
- ❑ Be attentive to manufacturer's expiration dates – don't re-invent reagent stability
- ❑ Storage of diluted standards should be validated
 - Storage containers
 - Storage temperature
 - Length of storage

Instrument Control

9

- Instruments must be properly qualified
 - Water baths/heat blocks mapped for heat distribution
 - Incubating plate/tube readers checked for temperature distribution and control, optics, data collection/transfer, data management, calculations
 - Some manufacturers calculate averages off of raw onset times, others off of test results
 - Pipettors must be calibrated. For example, for a 10 μ L spike, if the pipettor is off by 20%, there could be a significant impact on reported spike recovery

Test Method Control

10

- Suitability testing (inhibition/enhancement)
 - Per compendia (BET chapter is harmonized)
 - Understand your product – many conditions can affect the aggregation of endotoxin in solution
 - Lack of spike recovery may require a change in lysate vendor and/or test method
 - Lack of spike recovery may require treatment of the sample (heat, ultrafiltration, etc)
 - Keep abreast of dosing or administration changes as reflected in the package insert that could affect the endotoxin limit
 - If the endotoxin limit and MVD change as the result of the PI change, the method may need re-qualification

Standard Curve Control

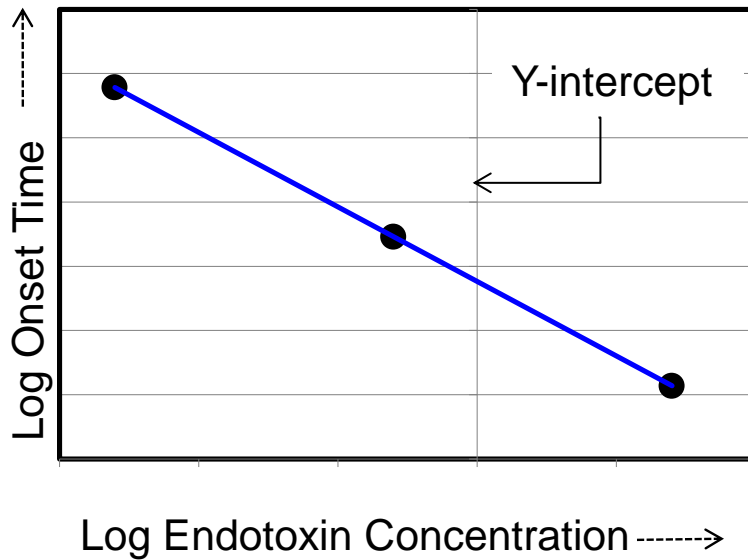
11

- Attributes that affect accuracy
 - Slope
 - Y-intercept
 - Linearity
 - % cv

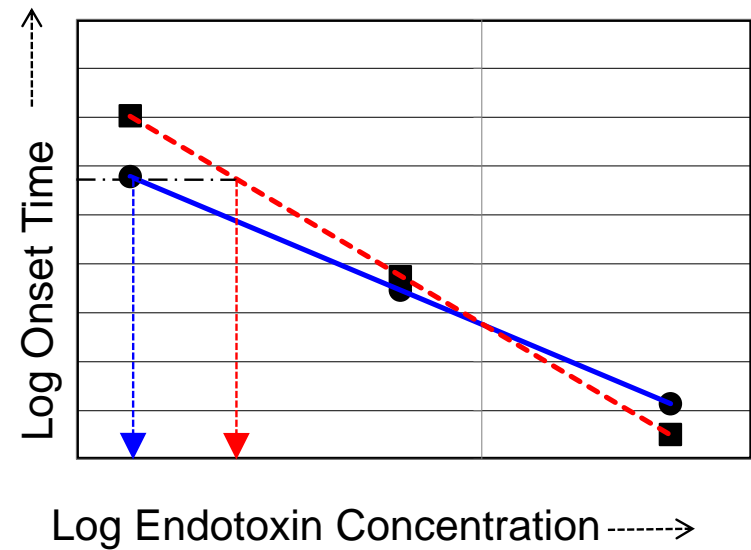
Standard Curve Control

12

“Perfect” curve



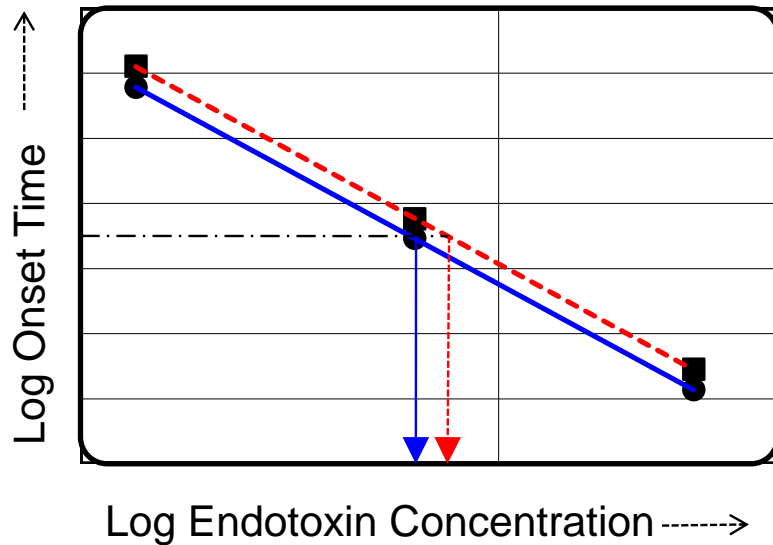
Change in slope



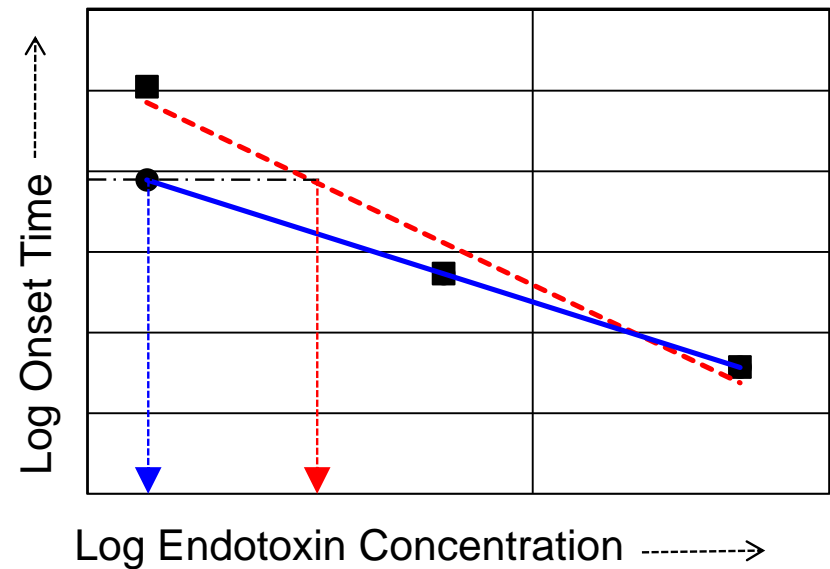
Standard Curve Control

13

1% change in y-intercept



$|r| = 0.982$



Data Interpretation

14

- What is an invalid test and what is OOS?
 - What happens on a quantitative test where one replicate is over the limit and one is under the limit and the average is OK?
- Trend invalid tests to identify issues in the laboratory CAPA
 - By product
 - By invalidity
 - By analyst
- Trend test results to identify adverse trends in manufacturing

Acme Pharmaceuticals 2012 BET Excursions and Invalid Tests

Date	Problem	Lysate Lot	Analyst	Product	correlation
1/1/12	invalid spike	XYZ123	BA	WFI port 2	0.991
1/15/12	OOS	XYZ123	KS	holding tank	0.998
2/12/12	void curve	XYZ123	BA	n/a	0.979
2/14/12	hot well	XYZ123	CC	NSA	0.999
2/22/12	invalid spike	XYZ123	BA	lipid emulsion	0.995
3/17/12	OOS	XYZ123	CC	ampicillin	0.999
4/1/12	invalid spike	XYZ123	BA	NSA	0.986
4/5/12	negative cont	XYZ123	MG	NaCl	0.996
5/1/12	hot well	XYZ123	CC	WFI port 2	0.999
5/9/12	OOS	XYZ123	BA	holding tank	0.997
5/29/12	invalid spike	ABC234	BA	Ringer's	0.993
6/8/12	invalid spike	ABC234	MG	lipid emulsion	0.999
6/14/12	hot well	ABC234	CC	recrystallized	0.986
7/5/12	OOS	ABC234	MG	holding tank	0.999
8/30/12	missed well	ABC234	KS	ampicillin	0.994
9/3/12	void curve	ABC234	BA	Ringer's	0.976
9/13/12	negative cont	ABC234	KS	WFI port 2	0.998
10/7/12	hot well	ABC234	CC	NaCl	0.999
10/31/12	invalid spike	ABC234	BA	holding tank	0.991
11/5/12	invalid spike	ABC234	CC	lipid emulsion	1.000
11/25/12	OOS	ABC234	KS	holding tank	0.998
12/15/12	negative cont	ABC234	CC	NSA	0.997
12/31/12	void curve	ABC234	BA	NaCl	0.975

Analysis: Acme Pharmaceuticals 2012 BET Excursions and Invalid Tests. Sorted by problem and analyst

Date	Problem	Lysate Lot	Analyst	Product	Correlation
2/14/12	hot well	XYZ123	CC	NSA	0.999
5/1/12	hot well	XYZ123	CC	WFI port 2	0.999
6/14/12	hot well	ABC234	CC	recrystallized	0.986
10/7/12	hot well	ABC234	CC	NaCl	0.999
1/1/12	invalid spike	XYZ123	BA	WFI port 2	0.991
2/22/12	invalid spike	XYZ123	BA	lipid emulsion	0.995
4/1/12	invalid spike	XYZ123	BA	NSA	0.986
5/29/12	invalid spike	ABC234	BA	Ringer's	0.993
10/31/12	invalid spike	ABC234	BA	holding tank	0.991
6/8/12	invalid spike	ABC234	MG	lipid emulsion	0.999
11/5/12	invalid spike	ABC234	CC	lipid emulsion	1.000
8/30/12	missed well	ABC234	KS	ampicillin	0.994
4/5/12	negative cont	XYZ123	MG	NaCl	0.996
9/13/12	negative cont	ABC234	KS	WFI port 2	0.998
12/15/12	negative cont	ABC234	CC	NSA	0.997
1/15/12	OOS	XYZ123	KS	holding tank	0.998
3/17/12	OOS	XYZ123	CC	ampicillin	0.999
5/9/12	OOS	XYZ123	BA	holding tank	0.997
7/5/12	OOS	ABC234	MG	holding tank	0.999
11/25/12	OOS	ABC234	KS	holding tank	0.998
2/12/12	void curve	XYZ123	BA	n/a	0.979
9/3/12	void curve	ABC234	BA	Ringer's	0.976
12/31/12	void curve	ABC234	BA	NaCl	0.975

Thank You!!!!!!

17

QUESTIONS?