



Conducting Effective OOS or OOT Investigations for Unexpected Results from the BET Assay

Dealing with Unexpected Results in Routine Testing

- In a well controlled aseptic production process, the likelihood of product adulteration by endotoxin should be minimal
- This means that the vast majority of results generated in the BET should be either below the Endotoxin Release Limit (ERL) for the product or undetectable i.e. below the detection limit of the assay
- However, unexpected results can and do occur with some frequency in QC laboratories, worldwide



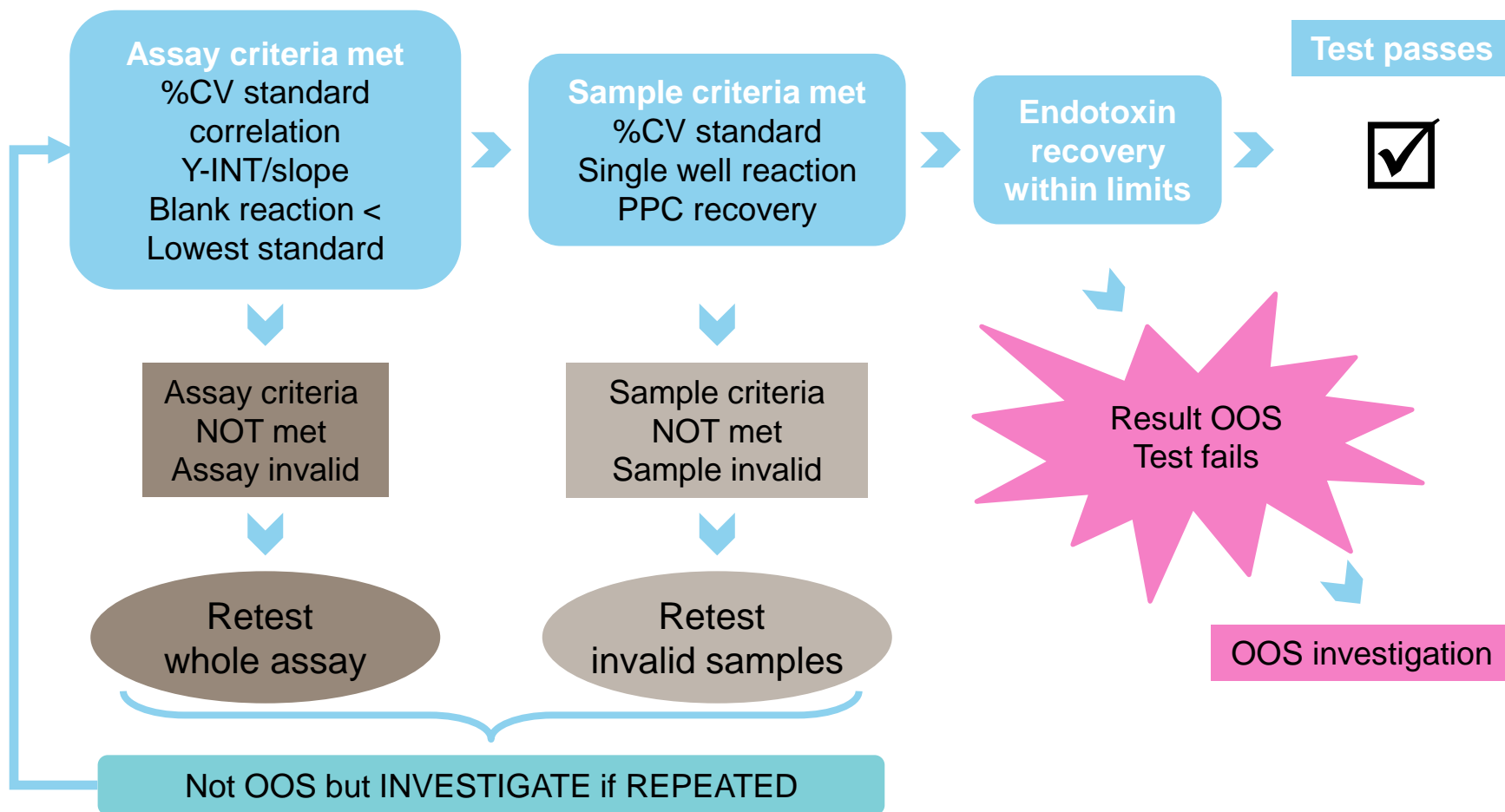
FDA Guidelines on OOS & OOT Results

- OOS (Out of Specification)
 - “All suspect results that fall outside the specification or acceptance criteria established in new drug applications, official compendia, or by the manufacturer”
- OOT (Out of Tolerance)
 - Consideration should equally be applied to results which are atypical and could constitute an increased risk
- Source
 - *FDA Guidance for Industry. Investigating Out of Specification (OOS) Test Results for Pharmaceutical Production.*
 - <http://www.fda.gov/cder/guidance/index.htm>

Initial Assessment

- There should be an objective and timely assessment to:
 - Establish likelihood of laboratory error
 - OR
 - Confirm there is an indication of manufacturing process / product problems
 - Assess significance of the result in overall QA program
- The process for making the decision on these key questions must be documented in an approved SOP

BET Assay Assessment Flow Chart



If an OOS Investigation Is Merited

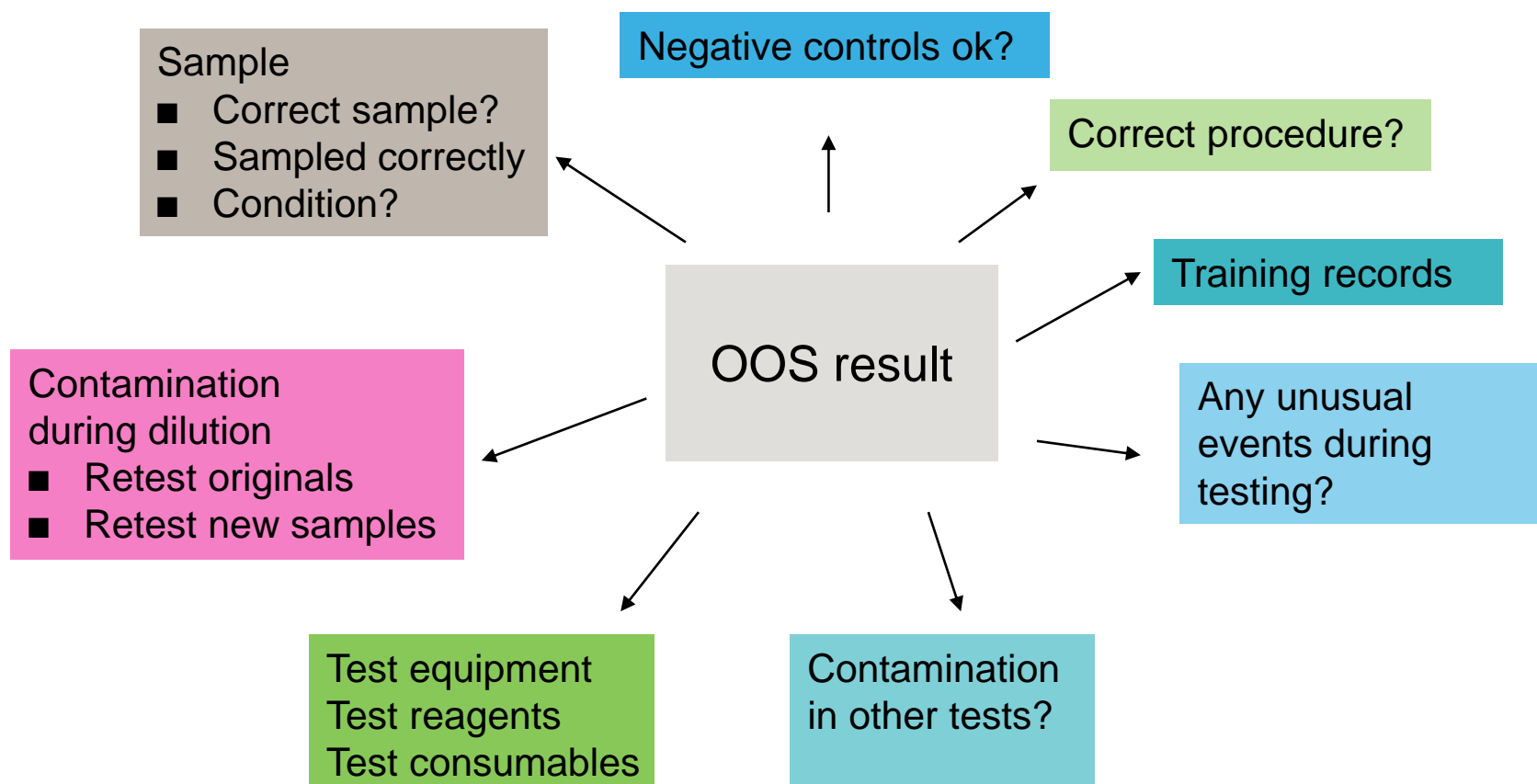
- Report suspect result to QA/Regulatory for assessment
- Notify production of need to investigate process
- An investigation team should be formed
- Start a documented investigation process, as soon as possible
- Make preliminary assessment of effects
- Inform customer services / production of potential consequences

Can the Result Be Attributed to Laboratory Error ?

- Supporting historical data with trend analysis will help
 - Justify the decision making process
 - Suggest Corrective And Preventative Actions (CAPA)
- Working with paper-based reports or excel spreadsheets can be cumbersome
- As kinetic BET assays generate large volumes of data, a database approach is recommended with Structured Query Language capability



Laboratory Error – Potential Causes



A Logical Approach Can Focus the Investigation

- When investigating potential laboratory error, use logic to narrow the field of investigation
- I usually classify issues with kinetic assays into one of three types
 - Likely to cause individual well variations, i.e. one well in error in a pair
 - Likely to cause multiple replicate errors, i.e. duplicate wells are in error
 - Likely to cause multiple well variations (all or several wells on the same plate in error)
- Similarly, you can identify factors that are likely to only affect the standard curve or only likely to affect sample results.
- Doing this will narrow the field of search, when asking questions of the database.

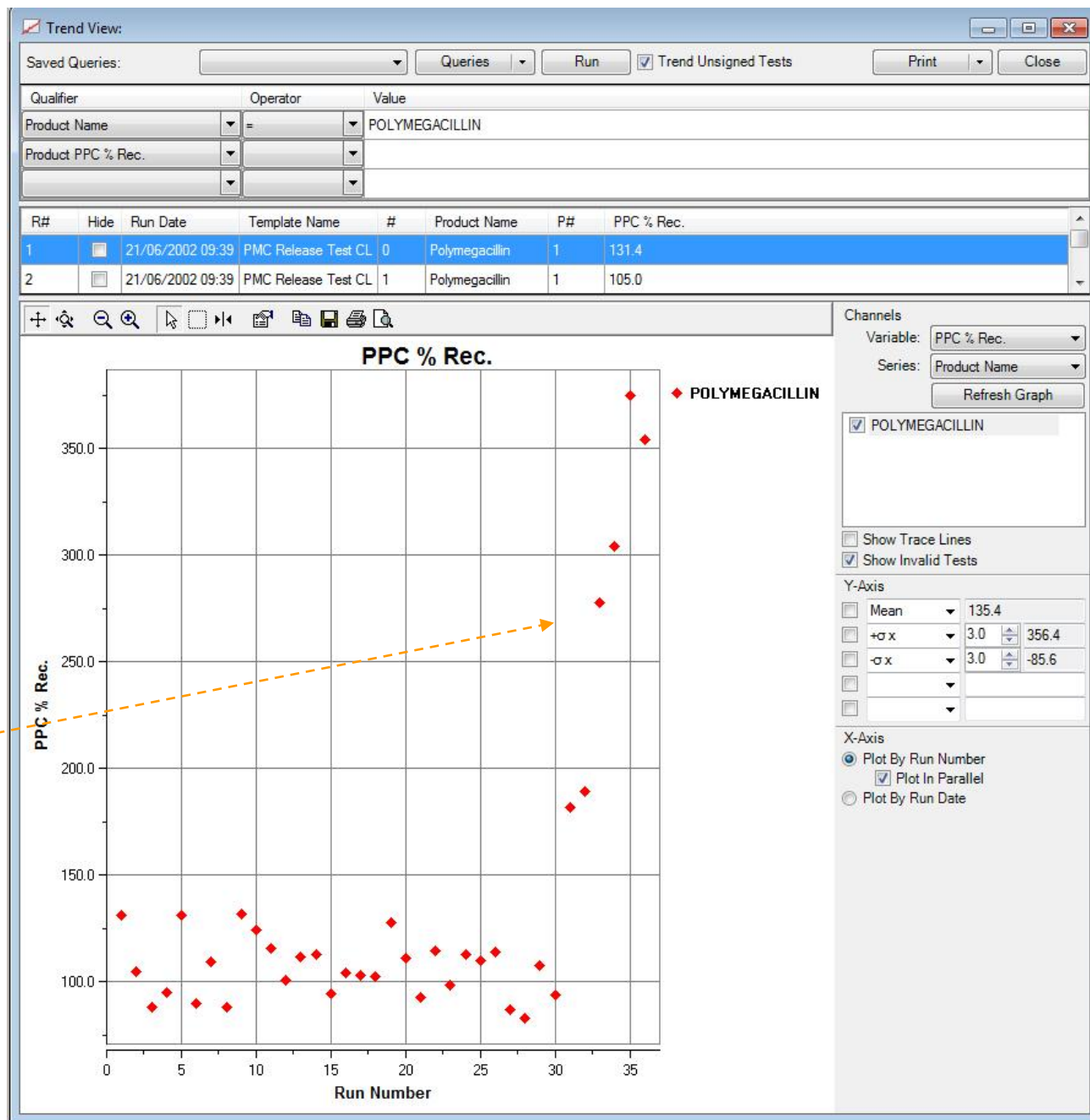
PPC Out-Of-Trend – Situation

- The following example illustrates the ease and efficiency of investigating OOS or OOT results using trend analysis:
 - Company 'XYZ' makes an antibiotic 'Polymegacillin'
 - PPC recovery percentages are normally consistent but recent assays show a sharp upward trend, with OOS & OOT results
 - Endotoxin results for the unspiked samples appear normal and follow previous trend

Potential Sources of Error in PPC Recovery

- What could be the possible cause(s)?
 - Technician Error?
 - Glucans?
 - Sample contamination?
 - Accessories?
 - Change in product formulation?
- Plan of action?
 - Data Analysis & Trending using the built-in Trend Analysis module of the WinKQCL™ Software

Query qualifiers used



Query qualifiers used
5 analysts involved

Trend View:

Saved Queries:

Queries

Run

☒ Trend Unsigned Tests

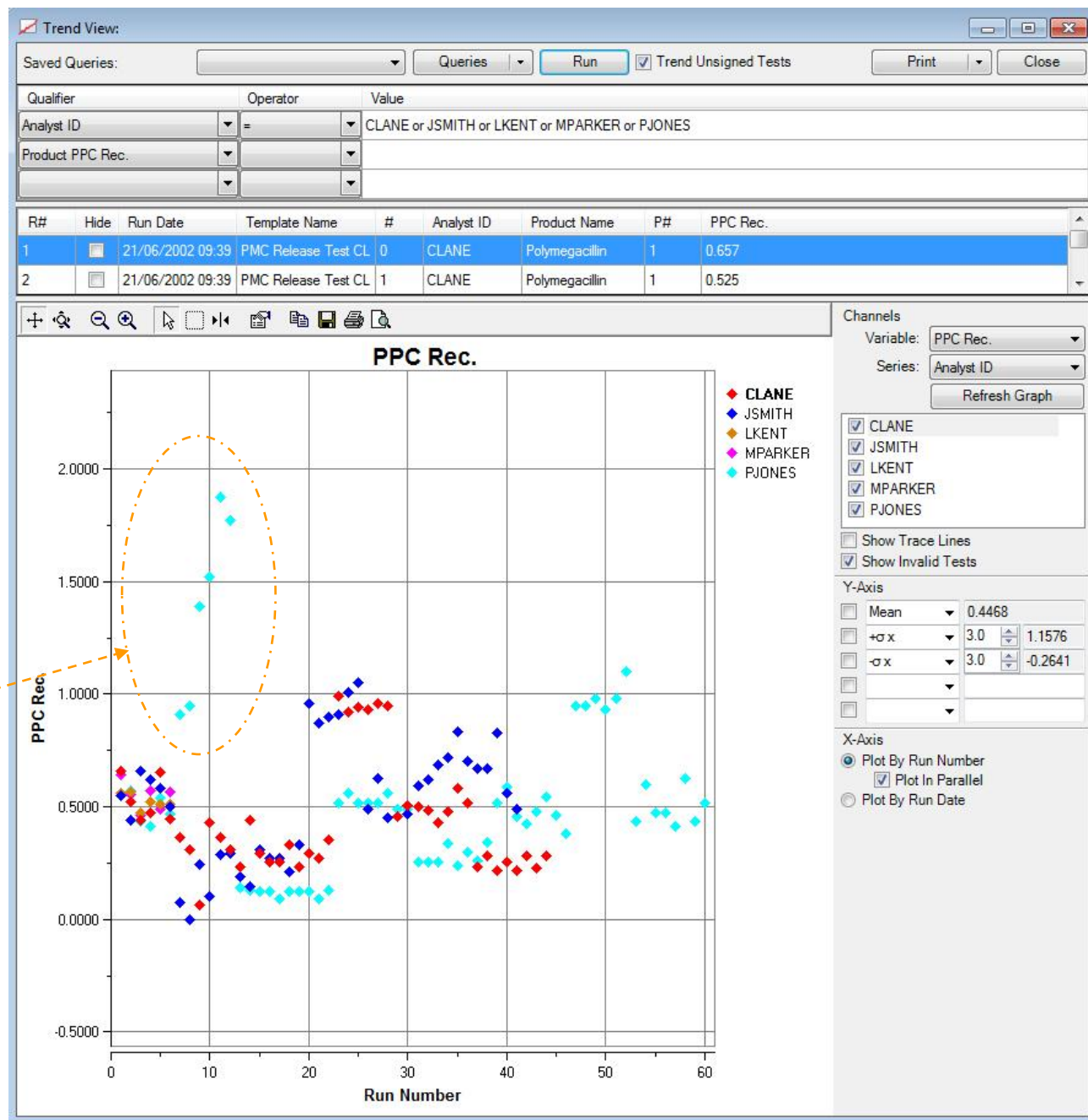
Print

Close

| Qualifier | Operator | Value |
|--------------|----------|----------------|
| Analyst ID | | |
| Product Name | = | POLYMEGACILLIN |
| | | |

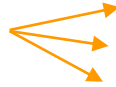
| R# | Run Date | Template Name | # | Analyst ID | Product Name | P# |
|----|------------------|----------------------|---|------------|----------------|----|
| 1 | 21/06/2002 09:39 | PMC Release Test CL | 0 | CLANE | Polymegacillin | 1 |
| 2 | 21/06/2002 09:39 | PMC Release Test CL | 1 | CLANE | Polymegacillin | 1 |
| 3 | 21/06/2002 09:39 | PMC Release Test CL | 2 | CLANE | Polymegacillin | 1 |
| 4 | 21/06/2002 09:39 | PMC Release Test CL | 3 | CLANE | Polymegacillin | 1 |
| 5 | 21/06/2002 09:39 | PMC Release Test CL | 4 | CLANE | Polymegacillin | 1 |
| 6 | 21/06/2002 09:39 | PMC Release Test CL | 5 | CLANE | Polymegacillin | 1 |
| 7 | 21/06/2002 10:27 | PMC Release Test JS | 0 | JSMITH | Polymegacillin | 1 |
| 8 | 21/06/2002 10:27 | PMC Release Test JS | 1 | JSMITH | Polymegacillin | 1 |
| 9 | 21/06/2002 10:27 | PMC Release Test JS | 2 | JSMITH | Polymegacillin | 1 |
| 10 | 21/06/2002 10:27 | PMC Release Test JS | 3 | JSMITH | Polymegacillin | 1 |
| 11 | 21/06/2002 10:27 | PMC Release Test JS | 4 | JSMITH | Polymegacillin | 1 |
| 12 | 21/06/2002 10:27 | PMC Release Test JS | 5 | JSMITH | Polymegacillin | 1 |
| 13 | 21/06/2002 11:10 | PMC Release Test LK | 0 | LKENT | Polymegacillin | 1 |
| 14 | 21/06/2002 11:10 | PMC Release Test LK | 1 | LKENT | Polymegacillin | 1 |
| 15 | 21/06/2002 11:10 | PMC Release Test LK | 2 | LKENT | Polymegacillin | 1 |
| 16 | 21/06/2002 11:10 | PMC Release Test LK | 3 | LKENT | Polymegacillin | 1 |
| 17 | 21/06/2002 11:10 | PMC Release Test LK | 4 | LKENT | Polymegacillin | 1 |
| 18 | 21/06/2002 11:10 | PMC Release Test LK | 5 | LKENT | Polymegacillin | 1 |
| 19 | 21/06/2002 13:46 | PMC Release Test ... | 0 | MPARKER | Polymegacillin | 1 |
| 20 | 21/06/2002 13:46 | PMC Release Test ... | 1 | MPARKER | Polymegacillin | 1 |
| 21 | 21/06/2002 13:46 | PMC Release Test ... | 2 | MPARKER | Polymegacillin | 1 |
| 22 | 21/06/2002 13:46 | PMC Release Test ... | 3 | MPARKER | Polymegacillin | 1 |
| 23 | 21/06/2002 13:46 | PMC Release Test ... | 4 | MPARKER | Polymegacillin | 1 |
| 24 | 21/06/2002 13:46 | PMC Release Test ... | 5 | MPARKER | Polymegacillin | 1 |
| 25 | 21/06/2002 14:28 | PMC Release Test PJ | 0 | PJONES | Polymegacillin | 1 |
| 26 | 21/06/2002 14:28 | PMC Release Test PJ | 1 | PJONES | Polymegacillin | 1 |
| 27 | 21/06/2002 14:28 | PMC Release Test PJ | 2 | PJONES | Polymegacillin | 1 |
| 28 | 21/06/2002 14:28 | PMC Release Test PJ | 3 | PJONES | Polymegacillin | 1 |
| 29 | 21/06/2002 14:28 | PMC Release Test PJ | 4 | PJONES | Polymegacillin | 1 |
| 30 | 21/06/2002 14:28 | PMC Release Test PJ | 5 | PJONES | Polymegacillin | 1 |

Query qualifiers used



Isolate items most
likely to be involved

e.g. P20 pipettor



Trend View:

Saved Queries:

Queries

Run

☒ Trend Unsigned Tests

Print

Close

| Qualifier | Operator | Value |
|----------------|----------|----------------|
| Product Name | = | POLYMEGACILLIN |
| Analyst ID | | |
| Accessory Name | | |
| | | |

| R# | Run Date | Template Name | # | Product Name | P# | Analyst ID | Accessory Name |
|----|------------------|---------------------|---|----------------|----|------------|---------------------|
| 1 | 21/06/2002 09:39 | PMC Release Test CL | 0 | Polymegacillin | 1 | CLANE | 13mm Dilution Tubes |
| 2 | 21/06/2002 09:39 | PMC Release Test CL | 1 | Polymegacillin | 1 | CLANE | 13mm Dilution Tubes |
| 3 | 21/06/2002 09:39 | PMC Release Test CL | 2 | Polymegacillin | 1 | CLANE | 13mm Dilution Tubes |
| 4 | 21/06/2002 09:39 | PMC Release Test CL | 3 | Polymegacillin | 1 | CLANE | 13mm Dilution Tubes |
| 5 | 21/06/2002 09:39 | PMC Release Test CL | 5 | Polymegacillin | 1 | CLANE | 13mm Dilution Tubes |
| 6 | 21/06/2002 09:39 | PMC Release Test CL | 4 | Polymegacillin | 1 | CLANE | 13mm Dilution Tubes |
| 7 | 21/06/2002 09:39 | PMC Release Test CL | 4 | Polymegacillin | 1 | CLANE | MP200 Multipettor |
| 8 | 21/06/2002 09:39 | PMC Release Test CL | 5 | Polymegacillin | 1 | CLANE | MP200 Multipettor |
| 9 | 21/06/2002 09:39 | PMC Release Test CL | 3 | Polymegacillin | 1 | CLANE | MP200 Multipettor |
| 10 | 21/06/2002 09:39 | PMC Release Test CL | 2 | Polymegacillin | 1 | CLANE | MP200 Multipettor |
| 11 | 21/06/2002 09:39 | PMC Release Test CL | 1 | Polymegacillin | 1 | CLANE | MP200 Multipettor |
| 12 | 21/06/2002 09:39 | PMC Release Test CL | 0 | Polymegacillin | 1 | CLANE | MP200 Multipettor |
| 13 | 21/06/2002 09:39 | PMC Release Test CL | 0 | Polymegacillin | 1 | CLANE | P20 Pipettor |
| 14 | 21/06/2002 09:39 | PMC Release Test CL | 1 | Polymegacillin | 1 | CLANE | P20 Pipettor |
| 15 | 21/06/2002 09:39 | PMC Release Test CL | 2 | Polymegacillin | 1 | CLANE | P20 Pipettor |
| 16 | 21/06/2002 09:39 | PMC Release Test CL | 3 | Polymegacillin | 1 | CLANE | P20 Pipettor |
| 17 | 21/06/2002 09:39 | PMC Release Test CL | 5 | Polymegacillin | 1 | CLANE | P20 Pipettor |
| 18 | 21/06/2002 09:39 | PMC Release Test CL | 4 | Polymegacillin | 1 | CLANE | P20 Pipettor |
| 19 | 21/06/2002 09:39 | PMC Release Test CL | 4 | Polymegacillin | 1 | CLANE | P200 Pipettor |
| 20 | 21/06/2002 09:39 | PMC Release Test CL | 5 | Polymegacillin | 1 | CLANE | P200 Pipettor |
| 21 | 21/06/2002 09:39 | PMC Release Test CL | 3 | Polymegacillin | 1 | CLANE | P200 Pipettor |
| 22 | 21/06/2002 09:39 | PMC Release Test CL | 2 | Polymegacillin | 1 | CLANE | P200 Pipettor |
| 23 | 21/06/2002 09:39 | PMC Release Test CL | 1 | Polymegacillin | 1 | CLANE | P200 Pipettor |
| 24 | 21/06/2002 09:39 | PMC Release Test CL | 0 | Polymegacillin | 1 | CLANE | P200 Pipettor |
| 25 | 21/06/2002 10:27 | PMC Release Test JS | 0 | Polymegacillin | 1 | JSMITH | 13mm Dilution Tubes |
| 26 | 21/06/2002 10:27 | PMC Release Test JS | 1 | Polymegacillin | 1 | JSMITH | 13mm Dilution Tubes |
| 27 | 21/06/2002 10:27 | PMC Release Test JS | 2 | Polymegacillin | 1 | JSMITH | 13mm Dilution Tubes |
| 28 | 21/06/2002 10:27 | PMC Release Test JS | 3 | Polymegacillin | 1 | JSMITH | 13mm Dilution Tubes |
| 29 | 21/06/2002 10:27 | PMC Release Test JS | 5 | Polymegacillin | 1 | JSMITH | 13mm Dilution Tubes |

Only two P20 units in active use

One unit PIP782 only used by technician PJONES

Trend View:

Saved Queries:

Queries

Run

☒ Trend Unsigned Tests

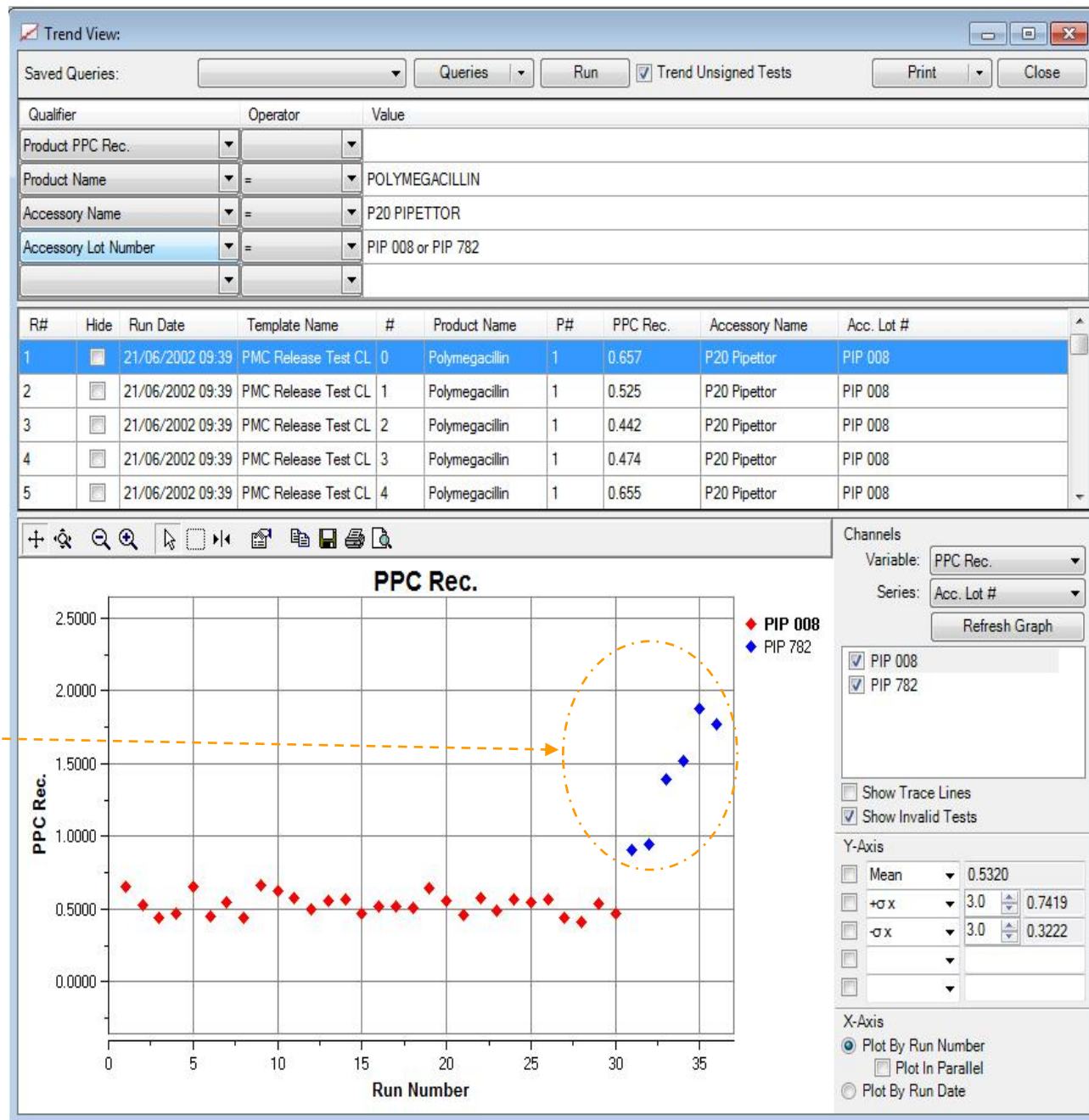
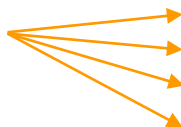
Print

Close

| Qualifier | Operator | Value |
|----------------------|----------|----------------|
| Product Name | = | POLYMEGACILLIN |
| Accessory Name | = | P20 PIPETTOR |
| Accessory Lot Number | | |
| Analyst ID | | |
| | | |

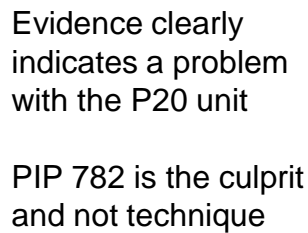
| R# | Run Date | Template Name | # | Product Name | P# | Accessory Name | Acc. Lot # | Analyst ID |
|----|------------------|----------------------|---|---------------|----|----------------|------------|------------|
| 19 | 21/06/2002 13:46 | PMC Release Test ... | 0 | Polmegacillin | 1 | P20 Pipettor | PIP 008 | MPARKER |
| 20 | 21/06/2002 13:46 | PMC Release Test ... | 1 | Polmegacillin | 1 | P20 Pipettor | PIP 008 | MPARKER |
| 21 | 21/06/2002 13:46 | PMC Release Test ... | 2 | Polmegacillin | 1 | P20 Pipettor | PIP 008 | MPARKER |
| 22 | 21/06/2002 13:46 | PMC Release Test ... | 3 | Polmegacillin | 1 | P20 Pipettor | PIP 008 | MPARKER |
| 23 | 21/06/2002 13:46 | PMC Release Test ... | 4 | Polmegacillin | 1 | P20 Pipettor | PIP 008 | MPARKER |
| 24 | 21/06/2002 13:46 | PMC Release Test ... | 5 | Polmegacillin | 1 | P20 Pipettor | PIP 008 | MPARKER |
| 25 | 21/06/2002 14:28 | PMC Release Test PJ | 0 | Polmegacillin | 1 | P20 Pipettor | PIP 008 | PJONES |
| 26 | 21/06/2002 14:28 | PMC Release Test PJ | 1 | Polmegacillin | 1 | P20 Pipettor | PIP 008 | PJONES |
| 27 | 21/06/2002 14:28 | PMC Release Test PJ | 2 | Polmegacillin | 1 | P20 Pipettor | PIP 008 | PJONES |
| 28 | 21/06/2002 14:28 | PMC Release Test PJ | 3 | Polmegacillin | 1 | P20 Pipettor | PIP 008 | PJONES |
| 29 | 21/06/2002 14:28 | PMC Release Test PJ | 4 | Polmegacillin | 1 | P20 Pipettor | PIP 008 | PJONES |
| 30 | 21/06/2002 14:28 | PMC Release Test PJ | 5 | Polmegacillin | 1 | P20 Pipettor | PIP 008 | PJONES |
| 31 | 21/06/2002 15:30 | PMC Release Test PJ | 0 | Polmegacillin | 1 | P20 Pipettor | PIP 782 | PJONES |
| 32 | 21/06/2002 15:30 | PMC Release Test PJ | 1 | Polmegacillin | 1 | P20 Pipettor | PIP 782 | PJONES |
| 33 | 21/06/2002 15:30 | PMC Release Test PJ | 2 | Polmegacillin | 1 | P20 Pipettor | PIP 782 | PJONES |
| 34 | 21/06/2002 15:30 | PMC Release Test PJ | 3 | Polmegacillin | 1 | P20 Pipettor | PIP 782 | PJONES |
| 35 | 21/06/2002 15:30 | PMC Release Test PJ | 4 | Polmegacillin | 1 | P20 Pipettor | PIP 782 | PJONES |
| 36 | 21/06/2002 15:30 | PMC Release Test PJ | 5 | Polmegacillin | 1 | P20 Pipettor | PIP 782 | PJONES |

Query qualifiers used



P20 with ID 782
is associated with
all high PPC values





Conclusions

- Use of a SQL-database trending program can help identify the cause of unexpected OOS or OOT results
- Such a database should be designed to automatically collect this data, as in the case in Lonza's WinKQCL™ Endotoxin Detection and Analysis Software
- The ability to cross-examine and trend data using query qualifiers helps to rapidly uncover the source(s) of error
- Use of such an integral database can save both time and money, whilst providing documentary evidence for decision processes

Thank You!