

Certificate of Analysis (example)

PRODUCT INFORMATION

Product: HEK-Blue™ Endotoxin Standard

Cat. code: rep-hbes

Lot number: O5E-XX-XX

Quantity: 50 EU

Storage temperature: -20°C

Recommended Retest Date: Month Year

QUALITY CONTROL

	Specifications	Results
Physico-chemical characteristics		
• <i>Appearance</i>	White lyophilized powder	Conform
• <i>Solubility (50 EU/ml in water for injection)</i>	Soluble	Conform
Endotoxin level		
• <i>Chromogenic LAL assay</i>	≥ 50 EU/mL	60 EU/mL
TLR2 contamination assay*		
• <i>TLR2 activity at 5 EU/ml</i>	Negative	Conform

* TLR activity tested on HEK-Blue™ TLR cells

Date: DD/MMM/YYYY

Reviewed by QA department:

Caution – Not fully tested. For research use only. Not for human use.