



## CERTIFICATE OF ANALYSIS

**Product Description:** PQ Low Control  
**Catalog Number/Size:** R92TE-Series  
**Lot Number:** Y0110082  
**Expiration Date:** 30 JUN 2023  
**Storage Temperature:** Room Temperature

Statement: PQ Low Control is verified to ensure that consistent signals are obtained across lots. Details on the quality control testing of PQ Low Control are provided in the product insert available at [www.mesoscale.com](http://www.mesoscale.com).

PQ Low Control was tested using SULFO-TAG™ Conjugated Biotinylated IgG (BTI) lot # R0010605 and MSD GOLD™ 96-well Streptavidin SECTOR Plate lot # Z0021486

Concentration of BTI	Mean ECL Spec and Tolerance Range	Intraplate CV or SD spec	Pass/Fail?
BTI with 1X	25565 - 47479	15 %	PASS
Dil-100 with 1X	0 - 50 counts	50 %	PASS

**Statement:** This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		21 AUG 2020