

## Product

Catalog Number: **CS2-0106-25mL**  
Description: **ViaStain™ AOPI Staining Solution**  
Lot Number: **200704-01-01**  
Appearance: **Orange**

Test	Result
AO Data Plot	Pass
PI Data Plot	Pass



**Matthew Bularzik, Quality Engineer**  
**07/15/2020**

Nexcelom Bioscience certifies that the above is an accurate record of the analysis of the subject lot. It is for research use only.