

Product Number:	102869
Lot Number:	503177
Expiration:	10-Apr-2022

Component	Part Number	Lot Number
Mouse IL-6 Sample Diluent	102104	110015
Mouse IL-6 Calibrator Diluent	102103	128012

Generic components- included with kit but not lot number specific		
Discovery Detector Diluent	101955	

Review/Approval

Shivani Goel QC Shivani Goel 21Oct2021
 Name Title Signature/Date