

Batch information

STD Mouse IFN gamma kit - Standard

Product reference : **62MIFNGCDA**

Product characteristics :

- Lot Number	03A
- Conditioning buffer	Phosphate 50 mM pH7.0 + BSA 1% + stabilizers
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2024.05.22

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	3 375.00 pg	N/A

PLEASE NOTE

Reconstitute with 0.375 mL H₂O.

The expiration date of batch 03A has been extended to 2024-05-22 after validation by QC.

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com

