

Batch information

STD HTRF Mouse IFN g kit - Std
Product reference : **62MIFNGCDA**

Product characteristics :

- Lot Number	05A
- Conditioning buffer	Phosphate 50 mM pH7.0 + BSA 1% + stabilizers
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2026.02.07

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	2 700.00 pg	N/A

PLEASE NOTE

Reconstitute with 0.300 mL H₂O.

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

