

## Batch information

**STD** HTRF Mouse IFN g kit - Std

Product reference : **62MIFNGCDA**

### Product characteristics :

- Lot Number	04A
- Conditioning buffer	Phosphate 50 mM pH7.0 + BSA 1% + stabilizers
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2025.01.26
- Reconstitute each vial with	0.3 mL distilled H2O

### 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](http://www.cisbio.com)*

