

## Batch information

### Mouse IFN $\beta$ kit- Standard

HTRF Mouse IFN beta kit - Std

Product reference :

**62MIFNBCDA**

### Product characteristics :

- Lot Number	01A
- Conditioning buffer	Phosphate 0.05M pH7 + 1% BSA + stabilizers
- Appearance	Lyophilized form
- Storage conditions	-20 °C
- Expiration date	2024.12.15
- Reconstitute each vial with	0.300000011920929 mL distilled H2O

### Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (pro)	1 800.00 pg	36 000

### PLEASE NOTE

The expiration date of batch 01A has been extended to 2024-12-15 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](http://www.cisbio.com)***

