

## Batch information

**Mouse STING WT bind kit - Std**

HTRF Mouse STING WT Bind. kit - Std

Product reference :

**64BDSTGMCD A****Product characteristics :**

- Lot Number	01A
- Conditioning buffer	Hepes buffer pH7 - BSA - preservatives
- Appearance	Frozen form
- Storage conditions	<-60 °C
- Expiration date	2024.08.23
- Volume per vial	50.00 µL
- Active Moiety Concentration (Pro)	2.00 µM

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Pro)	0.10 nmoles	N/A

**PLEASE NOTE**

The expiration date of batch 01A has been extended to 2024-08-23 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)***

