

## Batch information

**STD** Human TIMP1 kit - Standard

Product reference : **64TIMCDA**

### Product characteristics :

- Lot Number	01A
- Conditioning buffer	PO4 buffer pH7 - BSA - Preservatives
- Appearance	Frozen form
- Storage conditions	<-60 °C
- Expiration date	2023.05.15
- Volume per vial	150.00 µL
- Active Moiety Concentration (p)	250.00 ng/mL

### Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (p)	37.50 ng	N/A

### PLEASE NOTE

The expiration date of batch 01A has been extended to 2023-05-115 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)**

