

## Batch information

### STD

HTRF Human TIMP1 kit - Std

Product reference :

**64TIMCDA**

### Product characteristics :

|                                   |                                      |
|-----------------------------------|--------------------------------------|
| - Lot Number                      | 02A                                  |
| - Conditioning buffer             | PO4 buffer pH7 - BSA - Preservatives |
| - Appearance                      | Frozen form                          |
| - Storage conditions              | <-60 °C                              |
| - Expiration date                 | 2024.10.07                           |
| - 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

*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)*

