

Batch information

Tag-lite ligand 30

GLP1 receptor red agonist

Product reference :

L0030RED

Product characteristics :

| | |
|-----------------------|-------------|
| - Lot Number | 09A |
| - Conditioning buffer | DMSO |
| - Appearance | Frozen form |
| - Storage conditions | -20 °C |
| - Expiration date | 2023.11.01 |

Weight indications :

| | Quantity per vial | MW (Da) |
|---------------------|-------------------|---------|
| Active Moiety (Pro) | 0.46 nmoles | 4 999 |

PLEASE NOTE

The expiration date of batch 09A has been extended to 2023-11-01 after validation by CQ.

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com

