

Tag-lite ligand 30
Product reference :

GLP1 receptor red agonist
L0030RED

Product characteristics :

- Lot Number _____ 08A
- Conditioning buffer _____ DMSO
- Appearance _____ Frozen form
- Storage conditions _____ -20 °C
- Expiration date _____ 2022.05.01

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Pro)	0.47 nmoles	4 999

PLEASE NOTE

The expiration date of batch 08A has been extended to 2022-05-01 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