

Batch information

PY20-K

MAb PY20-Eu cryptate - 20,000 tests

Product reference :

61Y20KLB

Product characteristics :

- Lot Number	36A
- Packaging Unit	20 000 tests/vial
- Conditioning buffer	Hepes buffer 20mM pH7.4 + Stabilizers + 1%BSA
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2024.07.14
- Reconstitute each vial with	1 mL distilled H2O
- Active Moiety Concentration (Ab)	30.00 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	30.00 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (Ab)	1.50 ng
Number of tests/vial	20 000

*Actual amount per well should be optimized according to assay conditions

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