

## Batch information

**PY20-K**

MAb PY20-Eu cryptate - 5,000 tests

Product reference :

**61Y20KLA**

### Product characteristics :

- Lot Number	33A
- Packaging Unit	5 000 tests/vial
- Conditioning buffer	Hepes buffer 20mM pH7.4 + Stabilizers + 1%BSA
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2023.11.20
- Reconstitute each vial with	0.25 mL distilled H2O
- Active Moiety Concentration (Ab)	40.00 µg/mL

### Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	10.00 µg	150 000

### Recommended quantity per well\* :

	in 384 well low volume plate (20µL)
Active Moiety (Ab)	2.00 ng
Number of tests/vial	5 000

\*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](http://www.cisbio.com)**

