

**PY20-K**

MAb PY20-Eu cryptate - 5,000 tests

Product reference :

**61Y20KLA****Product characteristics :**

- Lot Number	32A
- 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	2023.10.29
- Reconstitute each vial with	0.25 mL distilled H2O
- Active Moiety Concentration (Ab)	50 µg/mL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Ab)	12.50 µg	150 000

**Recommended quantity per well\* :**

	in 384 well low volume plate (20µL)
Active Moiety (Ab)	2.50 ng
Number of tests/vial	5 000

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

**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)**