

Anti-IGE ETS07
Product reference :

Monoclonal antibody anti-IgE
MABETS07

Product characteristics :

- Lot Number_____ 01A
- Conditioning buffer_____ -
- Appearance_____ Lyophilized form
- Storage conditions_____ 2-8 °C
- Expiration date_____ 2028.07.03

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	100.00 µg	150 000

PLEASE NOTE

Reconstitute with 0.5 ml H₂O

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

