

Monoclonal antibody to Human IgA

Antibody:	MT57
Product code:	3860-3-250
Quantity:	250 µg
Immunogen:	Human IgA
Isotype:	Rat IgG2a
Specificity:	Human IgA. Cross-reacts with IgA from rhesus and cynomolgus macaques. Please visit www.mabtech.com for information on reactivity with other monkey species.
Concentration:	Supplied at 0.5 mg/ml in sterile filtered (0.2 µm) PBS with 0.02% sodium azide.
Purification:	Purified from <i>in vitro</i> cultures by protein G affinity chromatography.
Storage:	Store product at 4-8°C or frozen at -20°C or below. Avoid repeated freezing/thawing. The expiry date indicates how long unopened products, stored according to instructions, are recommended for use.
Applications:	For quantification of IgA in serum, plasma and cell culture supernatants using ELISA and for enumeration of IgA (total and antigen-specific) producing cells using ELISpot. Human IgA: In ELISA, MT57 is recommended as coating mAb in combination with ALP-conjugated detection mAb MT20 (product code 3860-9A). In ELISpot, MT57 is recommended as coating mAb in combination with biotinylated detection mAb MT20 (product code 3860-6). Monkey IgA: In ELISA and ELISpot MT57 is recommended as coating mAb in combination with anti-IgA-biotin (product code 3830-4).
Protocols:	Please contact Mabtech for ELISpot and ELISA protocols.



Note; for research use only.

Mabtech shall not be liable for the use or handling of the product or for consequential, special, indirect or incidental damages therefrom.

Mabtech AB (Head Office)
Sweden
Tel: +46 8 716 27 00
mabtech@mabtech.com

Mabtech, Inc.
USA
Tel: +1 513 871-4500
mabtech.usa@mabtech.com

Updated on 2020-03-26



Developed and manufactured by MABTECH AB, Sweden, whose quality management system complies with the standards ISO 9001:2015 & ISO 13485:2016.

www.mabtech.com