CTLA4 recombinant monoclonal antibody (060)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS974-0100

100µl

Manuals, SDS & CofA

View Online »

Handling & Storage

Handling Protected from light. Avoid freeze-thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Cytotoxic T-lymphocyte associated protein 4, CD152

Application ELISA

Clone 060

Formulation Liquid. 0.2 µm filtered solution in PBS. Preservative free.

Host Rabbit

Immunogen Recombinant human CTLA-4 / CD152 protein.

Isotype IgG

Purity Detail Protein A purified.

Recommendation Dilutions/Conditions ELISA (1:5'000-1:10'000)Suggested dilutions/conditions

may not be available for all applications. Optimal conditions

must be determined individually for each application.

Species Reactivity Human

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: September 17, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 info-

usa@enzolifesciences.com

European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com