## TIM-3 monoclonal antibody (6B6E2)

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

Order Online »

ENZ-ABS949-0100

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 T-cell immunoglobulin mucin receptor 3, HAVCR2,

Hepatitis A virus cellular receptor 2, TIMD3

**Application** ELISA

Clone 6B6E2

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

**Host** Mouse

**Immunogen** Recombinant human TIMD3 protein.

lsotype lgG2b

Purity Detail Protein A purified.

Recommendation Dilutions/Conditions ELISA (1:1'000-1:2'000) ELISA(Detection) (1:1'000-

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

Phone: 800.942.0430
infousa@enzolifesciences.com

SENZOLIFE SCIENCES
(PLS) AGE
(ELS) AG
phone: 441 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