

Data Sheet

Anti-OX40 Antagonist Antibody

Recombinant, Human Monoclonal
Anti-OX40

Catalog #: 72063-1

Lot #: 161005

Conc.: 0.79 mg/ml

Host Species | Isotype:

Human | Human IgG1

Formulated in: 8 mM Phosphate, pH 7.4,
110 mM NaCl, 2.2 mM KCl, and 20%
glycerol.

Stability: Stable for at least 12 months at
-80°C. *Protect from light. Avoid freeze/thaw
cycles. Store diluted antibody with carrier
protein (BSA 0.1 – 0.5%) at 4°C, stable for 2
weeks.*

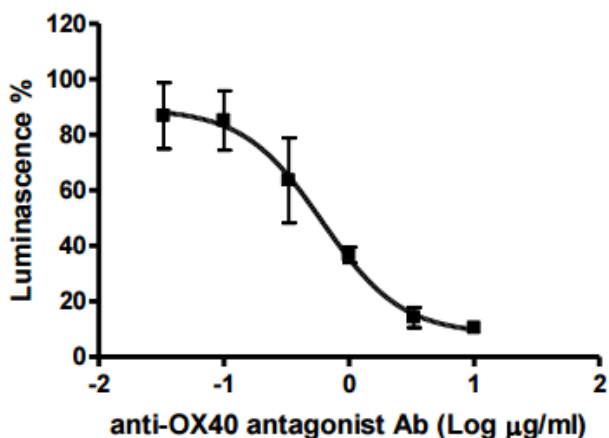
References:

1. Garrison, K., *et al.*, *Cancer Immunology*.
2012 April; **61(4)**: 511-521.
2. Bulliard, Y., *et al.*, *Immunology & Cell
Biology*. 2014 Apr; **92**: 475-480.

Description: Anti-OX40 recombinant human
antibody recognizing the OX40L binding region
of human OX40. This antibody has not been
tested for cross-reactivity with any other
species.

Purification: Protein A affinity chromato-
graphy from HEK293 supernatants.

Application: Useful as an antagonist to block
the signaling and binding of OX40 to its
ligand, OX40L in cellular and biochemical
assays, and for labeling cells expressing
OX40 for flow cytometry and
immunofluorescence microscopy.



Cells were treated with anti-OX40
antagonist antibody and 400 ng/ml of
OX40L simultaneously for 6 hours.

The IC₅₀ of anti-OX40 antagonist antibody
is 0.61 µg/ml

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com