

Data Sheet

Anti-OX40 Competitive Antibody

Recombinant, Human Monoclonal

Catalog #: 72558-2

Lot #: 160608

Conc.: 1.1 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.

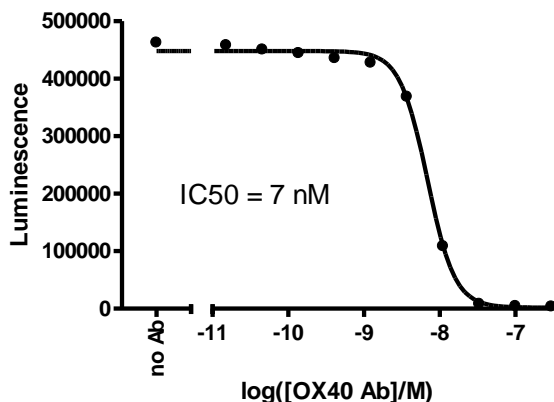
Description: Recombinant human Anti-OX40 competitive antibody recognizing the OX40 binding region. This antibody has not been tested for cross-reactivity with any other species.

Assay Conditions: Reaction done according to OX40[biotin]-OX40L assay kit (Catalog # 72045). Buffer is PBS + 0.1% BSA. Blocking buffer is Superblock + 0.05% Tween-20. Coat protein was added ON at 4 degrees. Inhibitor and then biotinylated ligand added for 2 hours at room temperature. Strep-HRP added for 1 hour at room temperature.

Purification: Protein A affinity chromatography from HEK293 supernatants.

Application: Useful for the inhibition of OX40:OX40L interactions.

OX40: OX40L Binding Assay



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