

Data Sheet

Anti-CD73 Antagonist Antibody

Recombinant, Human, Monoclonal
Catalog #: 79421-2
Lot #: 180205 **Conc.:** 1.06 mg/ml

Host Species | Isotype:

Human | IgG1

Formulated in: 50 mM Tris-HCl, pH 8.0
 138mM NaCl, 2.7 mM KCl.

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

Description: Anti-CD73 recombinant Human
 antagonist antibody. This antibody not been
 tested for cross-reactivity with other species.

Purification: Protein A affinity purification
 from HEK293 supernatants.

Application: Targeting CD73 has been show
 to prevent tumor recurrence and metastases
 in cancer patients. Anti-CD73 antibodies have
 been suggested as a new form of
 personalized cancer therapy.

References:

1. Deaglio, S. et al., J. Exp. Med. 2007;
 204(6): 1257-1265.
2. Thompson, L.F., et al., J. Exp. Med.
 2004; 200(11): 1395-1405.

Specific Activity

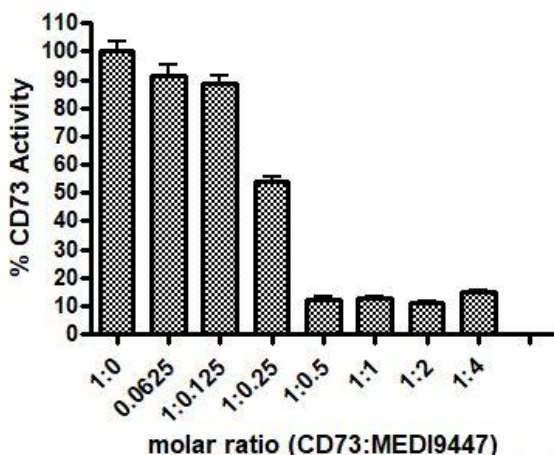


Fig.1 CD73 was incubated with anti-CD73 antibody in 1X CD73 reaction buffer (BPS #74000) at the indicated ratio for 1 hour at room temperature. Reaction was initiated by adding AMP (100 µM) and performed for 25 min at room temperature. CD73 activity was measured using Colorimetric Detection reagent (BPS #74001).

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