

Data Sheet

Anti-CD73 Antagonist Ab (clone #7F3)

Recombinant, Murine/Human Chimeric,
Monoclonal

Catalog #: 100231

Lot #: 181010 **Conc.:** 2.0 mg/ml

Host Species | Isotype:

Murine/Human Chimeric | IgG1

Formulated in: 40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl and 20% glycerol.

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-human CD73 recombinant Murine/Human chimeric 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 shown 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

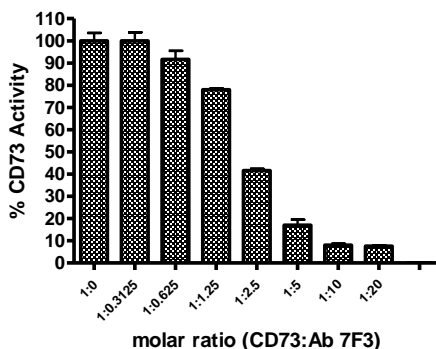


Fig.1 CD73 was incubated with anti-human CD73 antibody (7F3) in 1X CD73 reaction buffer (BPS #74000) at the indicated ratio for 1 hour at room temperature. Reaction was initiated by adding AMP (100 µM) and performed for 25 min at 37°C. CD73 activity was measured using Colorimetric Detection reagent (BPS #74001).

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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com