

Data Sheet

CD73, Avi-His-Tag, Biotin-Labeled (Mouse)

Mouse, Recombinant, C-terminal Avi-His-tag,
Biotin-Labeled

Catalog #: 79416

Lot #: 180315 **Conc.:** 0.34 mg/ml

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

Stability: At least 6 months at -80°C . Avoid
freeze/thaw cycles. Storing diluted protein is
not recommended, if necessary, use carrier
protein (BSA 0.1 – 0.5%).

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.

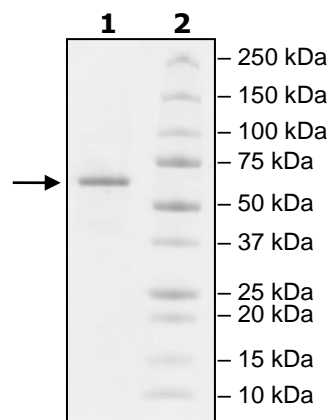
Description: Mouse CD73, also known as
5'-nucleotidase and Ecto-5'-nucleotidase,
GenBank Accession No. NM_011851, a.a.
29-549 with C-terminal Avi-His-tag,
expressed in a HEK293 cell expression
system. MW = 61 kDa. This protein runs at
a higher MW by SDS-PAGE due to
glycosylation. This protein is enzymatically
biotinylated using Avi-Tag™ technology.

Quality Assurance

**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
2µg CD73
Lane 2:
Protein Marker

MW: 61 kDa + glycans
Purity: ≥90%



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