

Data Sheet

CD73, Avi-His-Tag, Biotin Labeled

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

Catalog #: 79127

Lot #: 170817

Conc.: 0.36 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: Human CD73, also known as
5'-nucleotidase and Ecto-5'-nucleotidase,
GenBank Accession No. NM_002526, a.a.
27-547 with C-terminal Avi-His-tag,
expressed in a HEK293 cell expression
system. This protein is enzymatically
biotinylated using AviTag™ technology.
MW = 61 kDa.

Specific Activity: >20 nmol/min/μg

Assay Conditions: Experiment ran
according to CD73 colorimetric protocol (15
min incubation at 37°C). Phosphate
standard (from Enzo #BML-AK111) was
used at 100 μM serially diluted half-fold
down the plate. CD73 [B] used at 0.3 ng/μl
and serially diluted half-fold.

Application: Useful for the study of enzyme
kinetics, screening inhibitors, and selectivity
profiling.

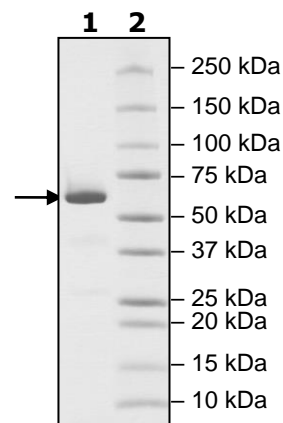
Quality Assurance

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

Lane 1:
2.5 μg CD73

Lane 2:
Protein Marker

MW: 61 kDa
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