

Data Sheet

CD73, Avi-His-Tag, Biotin Labeled

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

Catalog #: 79127

Lot #: 200122

Conc.: 0.31 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.

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.
Biotinylation is confirmed to be $\geq 90\%$. MW =
61 kDa.

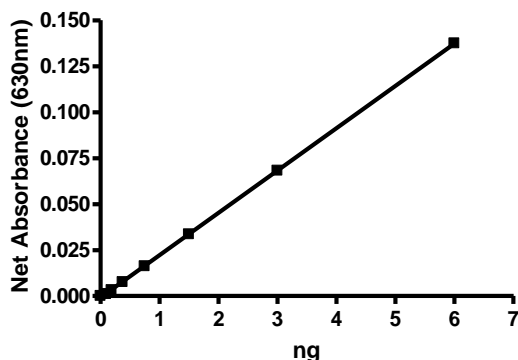
Specific Activity: >7000 pmol/min/μg

Assay Conditions: Assay was done in 50 μL
assay buffer containing 25 mM Tris, pH 7.4,
5 mM MgCl₂ and 100 μM AMP for 15 min at
37°C. Phosphate standard solution was
run side-by-side for product quantification.

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

Quality Assurance

Specific Activity:

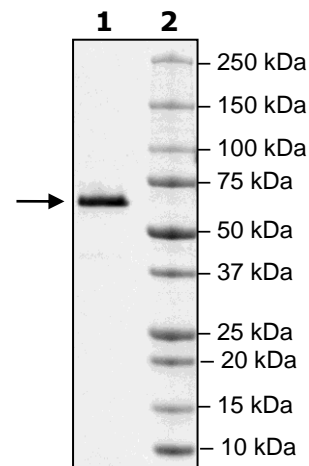


**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
2μg 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