

Data Sheet

BirA

Recombinant, N-terminal His-FLAG tag

Catalog #: 70032

Lot#: 130925-1 Conc.: 2.2 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, 0.04% Tween-
20, 20% glycerol, and 80 mM imidazole.

Stability: >6 months at -70°C . Avoid
freeze/thaw cycles.

References:

1. Li, Y., *et al.*, *Protein Expr Purif.* 2012
Mar; **82(1)**: 162-7.
2. Chakravarty, V., *et al.*, *J Bacteriol.*
2012 Mar; **194(5)**: 1113-26.

Description:

Biotin ligase BirA, GenBank Accession
No. AP012306, a.a. 2 – 321 (end) with
N-terminal His-FLAG tag, MW = 37
kDa, expressed in a baculovirus
infected Sf9 cell expression system.

Application:

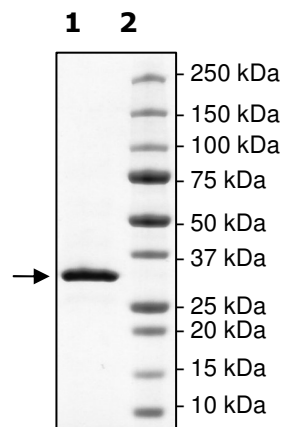
Useful for covalent *in vitro* biotin-
labeling of receptor proteins for
detection and purification assays.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
3.3 μg BirA
Lane 2:
Protein Marker

Purity: $\geq 90\%$
MW: 37 kDa



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



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence:

MHHHHHHDYKDDDDK KDNTVPLKLIALL ANGEFHSGEQLGETLGMSRAAINKHIQTL
RDWGVDFVFTVPGKGYSLPEPIQLLNAKQILGQLDGGSVAVLPVIDSTNQYLLDRIGEL
KSGDACIAEYQQAGRGRGRKWFSPFGANLYLSMFWRLEQQPAAIGLSLVIGIVMA
EVLRLGADKVRVKWPNDLYLQDRKLAGILVELTGKTGDAAQIVIGAGINMAMRRVE
ESVYNQGWITLQEAGINLDRNTLAAMLIRELRAALELFEQEGLAPYLSRWEKLDNFN
RPVKLIIGDKEIFGISRGIDKQGALLLEQDGIIPWMGGEISLRS AEK

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