

# BirA, His-FLAG-Tags (E. coli) (E. coli-derived) Recombinant

Catalog: 70030

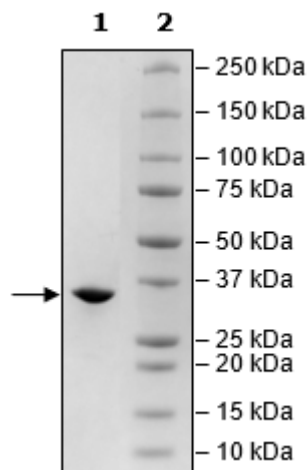
Lot: 220824

## Product Information

<b>Construct:</b>	BirA (His-FLAG-2-321(end))
<b>Concentration:</b>	1.13 mg/ml
<b>Species:</b>	<i>E. Coli</i>
<b>Formulated In:</b>	50 mM Tris-HCl, pH 8.0, 50 mM NaCl, 150 mM imidazole, 3 mM DTT, and 5% glycerol
<b>Expression System:</b>	<i>E. coli</i>
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	NP_418404
<b>MW:</b>	37 kDa
<b>Purity:</b>	≥90%
<b>Assay Conditions:</b>	BirA activity was confirmed using in vitro biotinylation of CBL-B by Western Blot. Lane 1: Standard; Lane 2: CBL-B, His-Avi-Tag (BPS Bioscience #80414); Lane 3: CBL-B, His-Avi-Tag (BPS Bioscience #80414) and BirA His-FLAG-Tags (E. coli) (BPS Bioscience 70030); Lane 4: CBL-B, His-Avi-Tag, Biotin-labeled (BPS Bioscience #80412). Protein loading was confirmed by Coomassie.
<b>Applications:</b>	Useful for covalently biotin-labeling of acceptor proteins for detection and purification assays

## Quality Control Data

### 4-20% SDS-PAGE Coomassie Staining



### BirA Activity

