

Data Sheet

KMO, His-FLAG-Tags

Human, Recombinant, C-terminal His-FLAG-tags

Catalog #: 11307

Lot #: 190225-1

Conc.: 0.46 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 100 µg/ml FLAG peptide, 0.2 mM TCEP and 20% glycerol.

Stability: At least 6 months at -80°C. *Avoid freeze/thaw cycles.*

References:

1. Zwilling, D., *et al.*, *Cell*, 2011; **6(10)**:863-874.
2. Giorgini, G., *et al.*, *Nature Genetics*. 2005; **37**: 526-531.

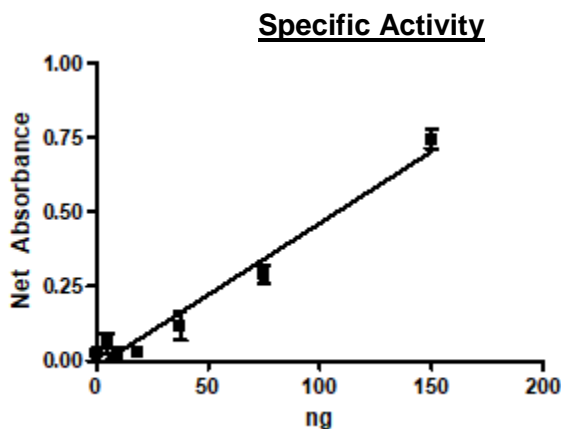
Description: Human KMO, also known as Kynurenine 3-Monooxygenase, Kynurenine 3-Hydroxylase, and EC 1.14.13.9, Genbank Accession No. NM_003679, a.a. 1-486 (end) with C-terminal His-FLAG-tag, expressed in a Baculovirus infected Sf9 cell expression system. MW = 58 kDa.

Specific Activity: 100 pmol/min/µg

Assay Conditions: KMO #11307 was tested with various amounts of enzyme. The reactions in KMO Assay Buffer (50 mM Sodium Phosphate, pH 7.5, 0.1% (w/v) Brij-35) contains 200 µM NADPH, 400 µM L-Kynurenine, and various concentrations of KMO. The reaction mixture incubated for 90 minutes while reading the UV absorption signal. For the negative control (blank), assay buffer was added instead of the enzyme. Absorption signals at 340 nm were measured using a Tecan Infinite M1000 plate reader.

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

Quality Assurance

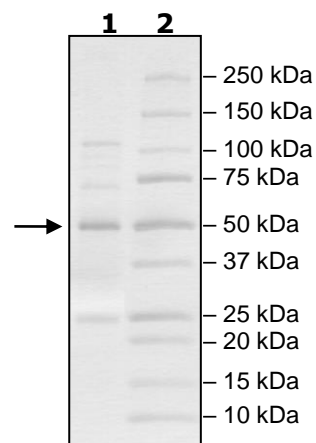


4-20% SDS-PAGE
Coomassie staining

Lane 1:
2 µg KMO

Lane 2:
Protein Marker

MW: 58 kDa
Purity: 60%



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