

Data Sheet

NOS2

Human, recombinant, N-terminal His-Flag tag

Catalog #: 70015

Lot#: 121018

Conc.: 0.15 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 80 ng/μl FLAG peptide, and 20% glycerol.

Stability: >6 months at -80 °C

References:

1. Hobbs, MR, *et al.*, *Lancet*. 2002 Nov 9;**360(9344)**:1468-75.
2. Okayama, H., *et al.*, *Int J Cancer*. 2012 May 23. doi: 10.1002/ijc.27644.

Description:

Human nitric oxide synthase 2 (NOS2) or inducible NOS (iNOS) (GenBank Accession No. NM_000625), amino acids 2 – 1153 (end) with N-terminal His-Flag tag, MW= 133 kDa, expressed in a baculovirus infected Sf9 cell expression system.

Application:

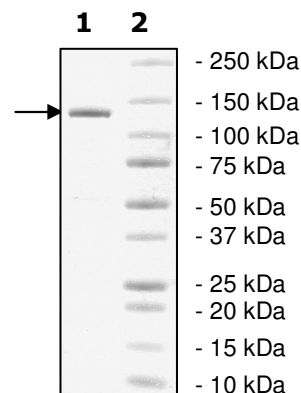
Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg NOS2
Lane 2:
Protein Marker

MW: 133 kDa
Purity: ≥97%



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 W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Protein Sequence:

MHHHHHHDYKDDDDK ACPWKFLFKTKFHQYAMNGEKDINNNVEKAPCATSSPVTQD
DLQYHNLSKQQNESPQPLVETGKKSPESLVKLDATPLSSPRHVRIKNWGSMTFQDT
LHHKAKGILTCRSKSLGSIMTPKSLTRGPRDKPTPPDELLPQAIEFVNQYYGSFKEA
KIEEHLARVEAVTKEIETTGTQYQLTGDELIFATKQAWRNAPRCIGRIQWSNLQVFDARS
CSTAREMFEHICRHVRYSTNNGNIRSAITVFPQRS DGKHDFRVWNAQLIRYAGYQMP
DGSIRGDPANVEFTQLCIDLGWKPKYGRFDVVPLVLQANGRDPELFEIPDLVLEVAM
EHPKYEFRELELKWYALPAVANMLLEVGGLEFPGCPFNWYMGTEIGVRDFCDVQ
RYNILEEVGRRMGLETHKLASLWKDQAVVEINI AVLHSFQKQNVTIMDHSAAESFMK
YMQNEYRSRGGCPADWIWL VPPMSG SITPVFHQEMLN YVLSPFY YQVEAWKTHVW
QDEKRRPKRREIPLKVLVKA VLFACMLMRKT MASRVRV TILFATETGKSEALAWDLG
ALFSCAFNPKVVCMDKYRLSCL EEEERLLL VVTSTFGNGDCPGNGEKLKKS LFM LKEL
NNKFRYAVFGLGSSMYPRFC AFAHDIDQKLSHLGASQLTPMGEGDEL SGQEDAFRS
WAVQTFKAACETFDVRGKQHIQIPKLYTSNVTWDPH HYRLVQDSQPLDLSKALSSMH
AKNVFTMRLKSRQNLQSPTSSRATILVELSCEDGQGLNYLPGEHLGVC PGNQPALVQ
GILERVVDGPTPHQTVRLEALDESGSYWVSDKRLPPCSLSQALTYFLDITTPPTQLLL
QKLAQVATEEPERQRLEALCQPSEYSKWKFTNSPTFLEVLEEFPSLRVSAGFLLSQL
PILKPRFYSISSSRDHTPTEIHLTVAVV TYHTRDGGGPLHHGVCSTWLNLSK PQDPVP
CFVRNASGFHLPEDPSHPCILIGPGTG IAPFRSFWQQRLHDSQHKGVRRGGRMTLVFG
CRRPDEDHIYQEEMLEMAQKGV LHAVHTAYSRLPGKPKVYVQDILRQQLASEVLRVL
HKEPGHLYVCGDVRMARDVAHTL KQLVAAKLKLN EEQVEDYFFQLKSQKRYHEDIF
GAVFPYEAKKDRVAVQPSSLEMSAL

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