

## Data Sheet

### NOS3

Human, recombinant, N-terminal His-Flag tag

**Catalog #:** 70018

**Lot#:** 121019      **Conc.:** 0.42 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 260 mM imidazole, and 20% glycerol.

**Stability:** >6 months at -80 °C

#### References:

1. Marsden, P.A., *et al.*, *FEBS Lett.* 1992 Aug 3;**307(3)**:287-93.
2. Liu, Y., Feng, Q., *Differentiation.* 2012 Jul;**84(1)**:54-61.

#### Description:

Human nitric oxide synthase 3 (NOS3), also known as endothelial NOS (eNOS), (GenBank Accession No. NM\_000603), amino acids 2 – 1203 (end) with N-terminal His-Flag tag, MW= 135 kDa, expressed in a baculovirus infected Sf9 cell expression system.

#### Application:

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

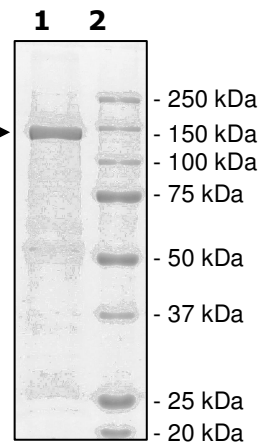
### Quality Assurance

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

**Lane 1:**  
1.6 µg NOS3 →

**Lane 2:**  
Protein Marker

**MW:** 135 kDa  
**Purity:** ≥71%



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](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://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](mailto:info@bpsbioscience.com)

**Protein Sequence:**

MHHHHHHDYKDDDDKGNLKSVAQEPGPPCGLGLGLGLGLGKQGPATPAPEPSRA  
PASLLPAPAHSPSSPLTQPPEGPKFPRVKNWEVGSITYDTLSAQAQQDGPCTPRR  
CLGSLVFPKRLQGRPSPGPPAPEQLLSQARDFINQYSSIKRSGSQAHEQRLQEVEA  
EVAATGTQYQLRESELVFGAKQAWRNAPRCVGRQWGLQVFDARDCRSAQEMFTYI  
CNHIKYATNRGNLRSAITVFPQRCPRGRDFRIWNSQLVRYAGYRQQDGSVRGDPANV  
EITELCIQHGWTGNGRFDVLPDLLQAPDDPELFLPPELVLEVPLEHPTLEWFAAL  
GLRWYALPAVSNMLLEIGGLEFPAAPFSGWYMSTEIGTRNLCDPHRYNILEDVAVCM  
DLDRTTSSLWKDKAAVEINVAVLHSYQLAKVTIVDHAATASFMKHLENEQKARGG  
CPADWAWIVPPISGSLTPVFHQEMVNYFLSPAIFYQDPWPKGSAKGTGTRKTKFK  
EVANAVKISASLMGTVMKRVKATILYGETGRAQSYAQLGRLFRKAFDPRVLCMD  
EYDVVSLEHETLVLVVTSTFGNGDPPENGESFAAALMEMSGPYNSSPRPEQHKS  
RFNSISCSPLVSSWRRKRKESNTDSAGALGTLRFVFLGSRAYPHFCFAFAVD  
TRLEELGGERLLQLGQDELGQEEAFRGWAQAQAFQAACETFCVGEDAKAAARDIF  
SPKRSWKRQRYRLSAQAEGQLLLPGLIHVHRRKMFQATIRSVENLQSSKSTRATILVR  
LDTGGQEGQLYQPGDHIGVCPNRPGLVEALLSRVEDPPAPTEPVAVEQLEKGS  
GPPGWVRDPRLLPCTLRQALTFFLDITSPSPQLLRLLSTLAEPPREQQELEALSQD  
PRRYEWWKWFRCPTLLEVLEQFSPVALPAPLLLTQLPLLQPRYYSVSSAPSTHPGEIH  
LTVAVLAYRTQDGLGPLHYGVCSTWLSQLKPGDPVPCFIRGAPSFRLPPDPSLPCILV  
GPGTGIAPFRGFWRERLHDIESKGLQPTMFLVFGCRCSQLDHLRDEVQNAQQRG  
VFGRLTAFSREPDNPKTYVQDILRTELAAEVHRLCLERGHMFCGDVTMATNVLQ  
TVQRILATEGDMELDEAGDVIGVLRDQQRVYHEDIFGLTLRTQEVTSRIRTSFSLQERQ  
LRGAVPWAFDPPGSDTNSP

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](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)