

Data Sheet

MUSK, His-Tag

Human, Recombinant, N-terminal His-tag
Catalog #: 40079
Lot#: 180416 **Conc.:** 0.40 mg/ml

Formulated in: 40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 3 mM DTT, 20% glycerol and 225 mM Imidazole.

Stability: At least 6 months at -80°C

References:

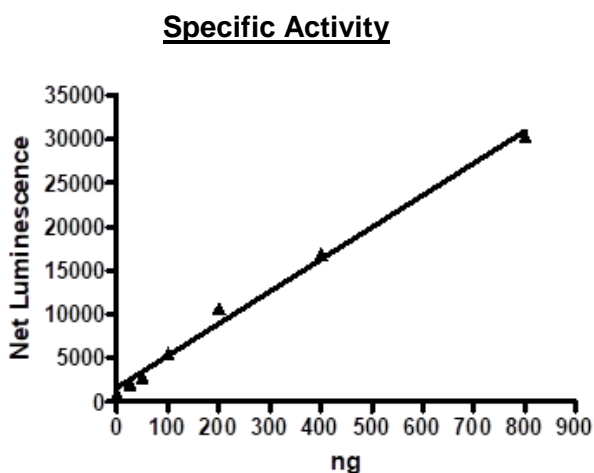
1. Zhou H. et al., J. Cell Biol. 146 (5), 1133-1146 (1999).
2. Evoli A. et al., Ann. N. Y. Acad. Sci. 1132, 76-83 (2008).

Description: Human MUSK, GenBank Accession No. NM_005592, a.a. 530-869 (end), with N-terminal His-tag, MW = 40 kDa, expressed in Baculovirus infected Sf9 cell expression system.

Assay Conditions: Assay was done using ADP-Glo (Promega) reagents with 100 μM ATP, 20 μM acetyl-CoA and 30 mM NaHCO₃ in 25 μL kinase assay buffer (30 mM HEPES pH=7.4, 2 mM MgCl₂, 1 mM DTT, and 2 mM Sodium citrate). Substrate used was MBP. Reaction was performed at 30°C for 40 min.

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

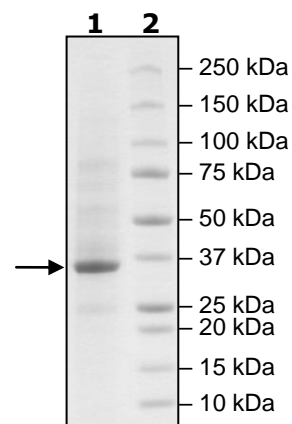
Quality Assurance



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

Lane 1:
2 μg MUSK
Lane 2:
Protein Marker

MW: 40 kDa
Purity: ≥66%



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 **1.858.259.5734**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com