

Product Information

Description:	Recombinant human NRF2 (Nuclear factor erythroid 2-related factor 2), full length, encompassing amino acids 3-605. This construct contains an N-terminal Flag tag. This recombinant protein was affinity purified.
Species:	Human
Construct:	NRF2 (FLAG-3-605(end))
Concentration:	1.26 mg/ml
Expression System:	<i>E. coli</i>
Purity:	≥90%
Format:	Aqueous buffer solution
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, and 3 mM DTT
MW:	69 kDa
Genbank Accession:	NM_006164.4
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.

Quality Control Data

4-20% SDS-Page Coomassie Staining

