

Product Information

Description:	Recombinant human NRF2 (Nuclear factor erythroid 2-related factor 2), full length, encompassing amino acids 3-605(end). This construct contains an N-terminal FLAG-tag. This recombinant protein was affinity purified.
Background:	Nuclear factor erythroid 2-related factor (Nrf2) is a transcription factor that coordinates antioxidant and anti-inflammatory responses by increasing the expression of antioxidant proteins through binding to the ARE (antioxidant response element) region of their promoters. Under basal conditions, Nrf2 is retained in the cytosol by cytoskeletal Keap1, a substrate recognition subunit of an E3 ligase complex which targets Nrf2 for degradation. Exposure to oxidative stress causes the release of Nrf2 from Keap1. Nrf2 translocates to the nucleus, where it binds to AREs and induces the expression of antioxidant and phase II proteins protecting the cell from oxidative damage. Both Nrf2 and Keap1 are attractive therapeutic targets.
Species:	Human
Construct:	NRF2 (FLAG-3-605(end))
Concentration:	0.28 mg/ml
Expression System:	<i>E. coli</i>
Purity:	86%
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.
Applications:	Useful for SDS-PAGE .

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

