

Recombinant IRF3 protein

Catalog No: 31544, 31944

Lot No: 31515001

Expressed In: Baculovirus

Quantity: 20 µg

Concentration: 0.3 µg/µl

Source: Human

Buffer Contents: Recombinant full length IRF3 protein is supplied at a concentration of 0.3 µg/µl in 25 mM HEPES pH 7.5, 300 mM NaCl, 5% glycerol, 0.04% Triton X-100, 0.2 mM TCEP

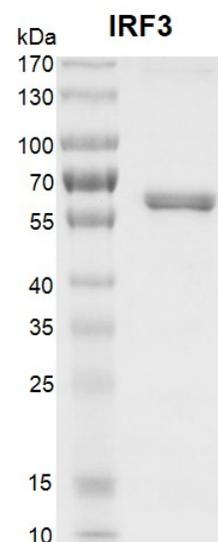
Background: IRF3 (Interferon regulatory factor 3) is a member of the IRF family of transcription factors that include IRF1, IRF6 and IRF7. IRF3 that has been phosphorylated at serine/threonine forms a complex with CREBBP and translocates to the nucleus to activate transcription of interferon-induced genes. IRF3 plays a critical role in the innate immune response against DNA and RNA viruses.

Protein Details: Recombinant human IRF3 was expressed in a baculovirus expression system as the full length protein (accession number NP_001562.1) with an N-terminal FLAG tag. The molecular weight of the protein is of 50.2 kDa. The IRF3 protein is >90% pure by SDS-PAGE.

Application Notes: This protein is useful for the study of protein-protein interaction, screening inhibitors, and selectivity profiling.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



Recombinant IRF3 protein gel.
IRF3 protein was run on a 10% SDS-PAGE gel and stained with Coomassie Blue.