

Recombinant SP140 (688-862) protein

Catalog No: 31448, 31848

Expressed In: *E. coli*

Quantity: 100, 1000 µg

Concentration: 1.5 µg/µl

Source: Human

Buffer Contents: Recombinant SP140 (688-862) protein was expressed in *E. coli* cells at a concentration of 1.5 µg/µl in 25 mM Tris pH 8.0, 500 mM NaCl, 0.04% Triton X-100, 20% glycerol.

Background: **SP140 nuclear body protein (SP140)** is a component of the nuclear body that may be involved in trafficking between the nucleus and the cytoplasm. SP140 contains one HSR domain, one PHD-type zinc finger, a bromodomain and one SAND domain. The bromodomain functions as a 'reader' of epigenetic histone marks and regulates chromatin structure and gene expression by linking associated proteins to the recognized acetylated nucleosomal targets. SP140 may be involved in the pathogenesis of acute promyelocytic leukemia and viral infection.

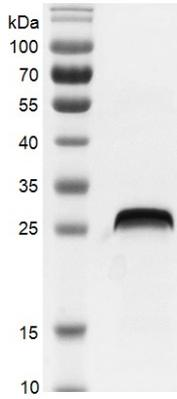
Protein Details: The peptide corresponding to amino acids 688-862 that contains the bromodomain sequences of SP140 (accession number NP_009168.4) was expressed in *E. coli* and contains an N-terminal His tag and C-terminal FLAG tag with an observed molecular weight of 26.98 kDa. The recombinant protein is ≥ 90% pure by SDS-PAGE.

Application Notes: Recombinant SP140 (688-862) is suitable for use in binding assays, inhibitor screening, 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.

SP140 (688-862)



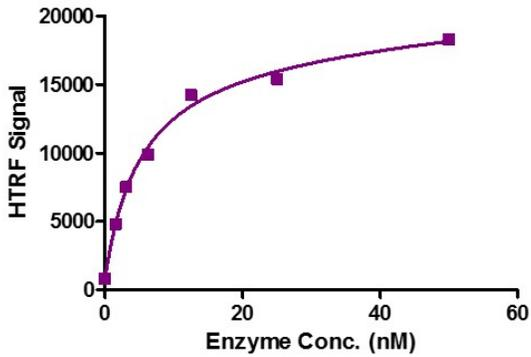
Recombinant SP140 (688-862) protein gel.

SP140 (688-862) protein was run on a 12.5% SDS-PAGE gel and stained with Coomassie Blue.

MW: 26.9 kDa

Purity: > 90%

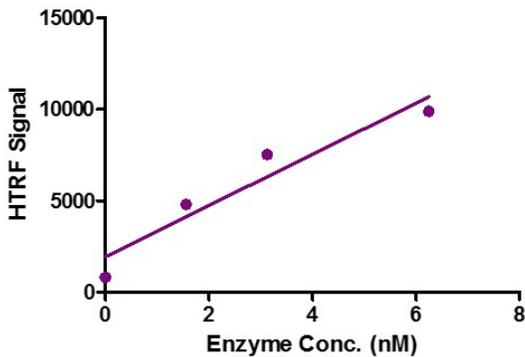
SP140 (688-862) Titration



HTRF Assay for Recombinant SP140 (688-862) protein activity.

3 μ M histone peptide H3K14ac was incubated with SP140 (688-862) in reaction buffer including 50 mM HEPES-NaOH pH 7.0, 0.1% BSA for 1 hour at room temperature. Anti-FLAG antibody was used to detect reaction products.

SP140 (688-862) Titration



HTRF Assay for Recombinant SP140 (688-862) protein activity.

3 μ M histone peptide H3K14ac was incubated with SP140 (688-862) in reaction buffer including 50 mM HEPES-NaOH pH 7.0, 0.1% BSA for 1 hour at room temperature. Anti-FLAG antibody was used to detect reaction products.