
FOSL2 Polyclonal Antibody

(Catalog # A71553)

Background

The Fos gene family consists of 4 members: FOS, FOSB, FOSL1, and FOSL2. These genes encode leucine zipper proteins that can dimerize with proteins of the JUN family, thereby forming the transcription factor complex AP-1. As such, the FOS proteins have been implicated as regulators of cell proliferation, differentiation, and transformation.

Description

FOSL2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Rat

Isotype

IgG

Uniprot ID

P15408

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 187-326 of human FOSL2 (NP_005244.1).

Storage

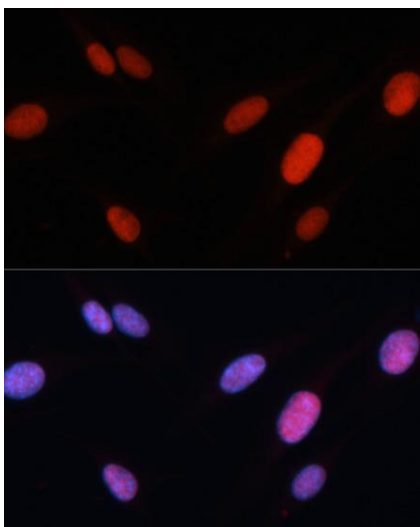
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

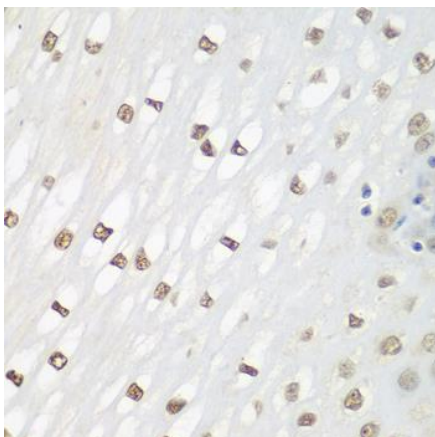
FOSL2; FRA2; fos-related antigen 2

Application

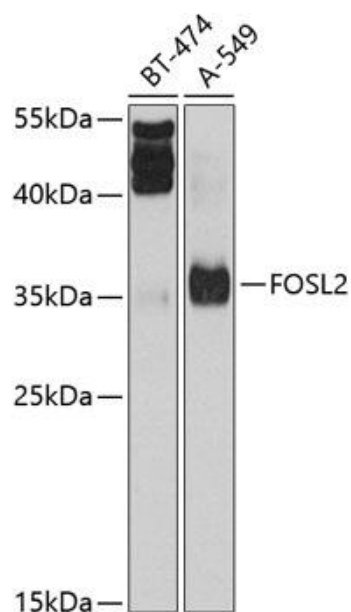
WB, IHC, IF; Recommended dilution: WB: 1:500-1:2000, IHC: 1:100-1:200, IF:1:50-1:200



Immunofluorescence analysis of U-2 OS cells using FOSL2 Polyclonal Antibody at dilution of 1:100. Blue: DAPI for nuclear staining.



Immunohistochemistry of paraffin-embedded human esophagus using FOSL2 Polyclonal Antibody at dilution of 1:100 (40x lens).



Western blot analysis of extracts of various cell lines, using FOSL2 Polyclonal Antibody at 1:1000 dilution. Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25ug per lane. Blocking buffer: 3% nonfat dry milk in TBST. Exposure time: 30s.