

---

## SGF29 Polyclonal Antibody, FITC Conjugated

(Catalog #A58448)

---

### Background

Involved in transcriptional regulation, through association with histone acetyltransferase (HAT) SAGA-type complexes like the TFTC-HAT, ATAC or STAGA complexes. Specifically recognizes and binds methylated 'Lys-4' of histone H3 (H3K4me), with a preference for trimethylated form (H3K4me3). In the SAGA-type complexes, required to recruit complexes to H3K4me. May be involved in MYC-mediated oncogenic transformation.

### Description

SGF29 Polyclonal Antibody, FITC Conjugated. FITC. Raised in: Rabbit.

### Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

### Specificity

Human

### Isotype

IgG

### Uniprot ID

Q96ES7

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human SAGA-associated factor 29 homolog protein (1-293AA)

### Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

SAGA-associated factor 29 homolog, CCDC101, SGF29, Coiled-coil domain-containing protein 101

### Application

ELISA, IHC; Recommended dilution: ELISA 1:100 - 1:500

---