

## Ski3 polyclonal antibody - Classic

**Cat. No. C15410011**

Type: Polyclonal	Specificity: Human
Size: 150 µg/50 µl	Isotype: NA
Concentration: 3.0 µg/µl	Host: Rabbit
Lot No.: 001	Purity: Protein G purified
Storage buffer: PBS containing 0.05% azide and 0.05% ProClin 300	Storage conditions: Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles.
Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.	

Last Data Sheet Update: November 5, 2019

### Description

Alternative names: **TTC37**

Polyclonal antibody raised in rabbit against human SKI3, using a recombinant protein.

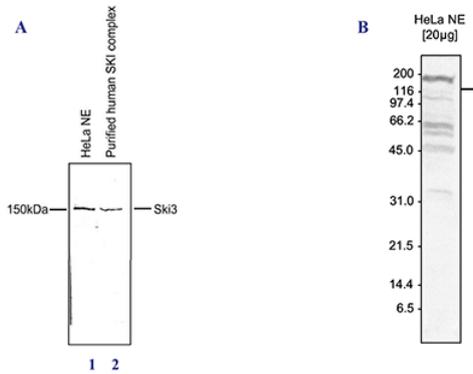
### Applications

Applications	Suggested dilution	References
Western Blotting	1:100	Fig 1
Immunofluorescence	1:100	Fig 2, Ref 1
ChIP	1 - 5 µg/ChIP	Ref 1

### Target Description

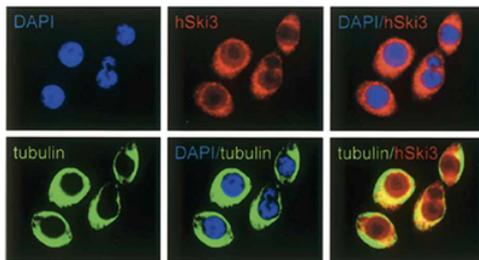
SKI3 is a component of the SKI complex which is composed of SKI2, SKI3 and SKI8. The SKI complex interacts with SKI7, making the link between the SKI complex and the exosome in order to perform mRNA degradation.

**Validation Data**



**Figure 1. Western blot analysis using the Diagenode antibody directed against SKI3**

Western blot was performed on HeLa nuclear extracts (lane 1) and on the purified human SKI complex (lane 2) (Figure 1A), or on nuclear extracts from HeLa cells (HeLa NE, 20 µg) (Figure 1B) using the Diagenode antibody against SKI3 (Cat. No. C15410011) diluted 1:100. The location of the protein is indicated on the right, the marker (in kDa) is shown on the left.



**Figure 2. Immunofluorescence with the Diagenode antibody directed against SKI3**

HeLa cells were stained with the Diagenode antibody against SKI3 (Cat. No. C15410011) (in red), with DAPI to label the DNA (in blue), and with an antibody against α-tubulin to label the cytoplasm (in green). SKI3 is detected in both the cytoplasm and the nucleus but appears to be excluded from the nucleolus.