

## A20

Cat. # DB209

### Background

A20, also known as TNFAIP3 (Tumor Necrosis Factor Alpha-Induced Protein 3), is a critical regulator of inflammation and immune responses in the human body. This protein acts as a dual-function enzyme with both deubiquitinase (DUB) and E3 ubiquitin ligase activities. A20's primary role is to dampen excessive immune responses by negatively regulating signaling pathways triggered by pro-inflammatory molecules, such as tumor necrosis factor-alpha (TNF- $\alpha$ ). It accomplishes this by removing ubiquitin chains from specific signaling molecules and adding inhibitory ubiquitin modifications to others. Dysregulation of A20 has also been associated with autoimmune diseases and inflammatory disorders, highlighting its importance in maintaining immune system balance.

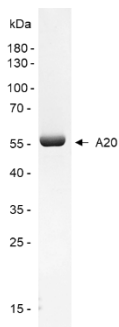
### Alternate names

TNF alpha-induced protein 3, OTU domain-containing protein 7C, Putative DNA-binding protein A20, Zinc finger protein A20

### Product Information

<b>Purity</b>	$\geq 95\%$ by SDS-PAGE
<b>Molecular Weight</b>	60 kDa
<b>Quantity</b>	25 $\mu$ g
<b>Physical State:</b>	Liquid
<b>Source</b>	Human Recombinant
<b>Tag</b>	His6
<b>Storage</b>	-80°C. Avoid repeated freeze/thaw cycles

### Product QC



**SDS-Page Analysis of purified A20.** Two  $\mu$ g of the enzyme was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.

### References

1. Momtazi, G., et al., Am J Physiol Lung Cell Mol Physiol., 2019. 316(3):L456-469.
2. Yu, M., et al., World J Pediatr., 2020. 16(6):575-584.

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2009 LifeSensors, Inc. All Rights Reserved