

A20/TNFAIP3

(Tumor necrosis factor alpha-induced protein 3, OTU domain-containing protein 7C, Putative DNA-binding protein A20, Zinc finger protein A20, OTUD7C, TNAP3)

- **CAT#:** A20/TNFAIP3
- **Enzyme:** Human A20/TNFAIP3
- **Substrate:** Fluorogenic substrate [Ubiquitin-AMC], 100 nM
- **Reference Compound IC50s:**

Compounds	IC50 (μ M)
UCHL1 I	41.8
Ub Aldehyde	5.64
UCHL1 I 2	65.8

- **Scientific Information:** P21580 (<http://www.expasy.org/uniprot/P21580>)

Reference Compound IC50 for A20/TNFAIP3

