

Ataxin-3/ATXN3

(Machado-Joseph disease protein 1, Spinocerebellar ataxia type 3 protein, ATX3, MJD, MJD1, SCA3)

- **CAT#:** Ataxin-3/ATXN3
- **Enzyme:** Human Ataxin-3/ATXN3
- **Substrate:** Fluorogenic substrate [Ubiquitin-AMC], 100 nM
- **Reference Compound IC50s:**

Compounds	IC50 (μ M)
UCHL1 I 2	153
UCHL3 I	298

- **Scientific Information:** P54252 (<http://www.expasy.org/uniprot/P54252>)

Reference Compound IC50 for Ataxin-3/ATXN3

