

## HDAC3/NCOR2

(Histone Deacetylase 3, Class I)

NCOR2: Nuclear receptor corepressor 2, CTG repeat protein 26, SMAP270)

- **CAT#:** HDAC3/NCOR2
- **Enzyme:** Human HDAC3/NCOR2
- **Substrate:** Fluorogenic peptide from p53 residues 379-382 (RHKK<sub>Ac</sub>), 50  $\mu$ M
- **Reference Compound IC50s:**

| Compounds      | IC50 (nM) |
|----------------|-----------|
| Trichostatin A | 11.3      |
| SAHA           | 132       |

- **Scientific Information:** HDAC3: O15379 (<http://www.expasy.org/uniprot/O15379>)  
NCOR2: Q9Y618 (<http://www.expasy.org/uniprot/Q9Y618>)

### Reference Compound IC50 for HDAC3

