

# HDAC Family Antibody Set

**CATALOG #:** K333-11-30

**AMOUNT:** 11 x 30 µg

**LOT #:** \_\_\_\_\_

**CONTENTS:**

Antibodies	Host	Species	Applications	Amount	Cap code	Part#
HDAC-1	Rabbit	hu/ms/rt	WB/IP/IHC	30 µg	Yellow	3601-30
HDAC-2	Rabbit	hu/ms/rt	WB/IP/IF	30 µg	Orange	3602-30
HDAC-3	Rabbit	hu/ms/rt	WB/IHC	30 µg	Red	3603-30
HDAC-4	Rabbit	hu/ms/rt	WB	30 µg	Orange	3604-30
HDAC-5	Rabbit	hu/ms/rt	WB/IP/IHC	30 µg	Blue	3605-30
HDAC-6	Rabbit	hu/ms/rt	WB/IHC	30 µg	Violet	3606-30
HDAC-7	Rabbit	hu/ms/rt	WB/IHC	30 µg	Red	3607-30
HDAC-8	Rabbit	hu/ms/rt	WB/IP/IHC	30 µg	Green	3608-30
HDAC-9	Rabbit	hu/ms/rt	WB/	30 µg	Blue	3609-30
HDAC-10	Rabbit	hu/ms/rt	WB	30 µg	Brown	3610-30
HDAC-11	Rabbit	hu/ms/rt	WB	30 µg	Yellow	3611-30

**FORMULATION:**

30 µg affinity purified rabbit anti-HDAC (1-11) polyclonal antibodies in phosphate buffered saline (PBS), pH 7.2, containing 30% glycerol, 0.5% BSA, 0.01% thimerosal.

**STORAGE CONDITIONS:**

Store at -20°C. For long-term storage, aliquot and freeze at -70°C. Avoid repeated freeze/thaw cycles.

**BACKGROUND:**

Histone deacetylase (HDAC) family proteins play an important role in modulating the eukaryotic chromatin structure.

**APPLICATION AND USAGE:**

The antibody can be used for Western blotting (0.5-4 µg/ml), Immunoprecipitation (10-20 µg/ml), and Immunohistochemistry (10-20 µg/ml). However, the optimal conditions should be determined individually. Blocking peptide for each antibody is available separately.

**NOTE:** For Research Use Only! Not to be used in human subjects.

**RELATED PRODUCTS:**

Apoptosis Detection Kits & Reagents

- Annexin V Kits & Bulk Reagents
- Caspase Assay Kits & Reagents
- Mitochondrial Apoptosis Kits & Reagents
- Nuclear Apoptosis Kits & Reagents
- Apoptosis Inducers and Set
- Apoptosis siRNA Vectors

Cell Fractionation System

- Mitochondria/Cytosol Fractionation Kit
- Nuclear/Cytosol Fractionation Kit
- Membrane Protein Extraction Kit
- Cytosol/Particulate Rapid Separation Kit
- Mammalian Cell Extraction Kit
- FractionPREP Fractionation System

Cell Proliferation & Senescence

- Quick Cell Proliferation Assay Kit
- Senescence Detection Kit
- High Throughput Apoptosis/Cell Viability Assay Kits
- LDH-Cytotoxicity Assay Kit
- Bioluminescence Cytotoxicity Assay Kit
- Live/Dead Cell Staining Kit

Cell Damage & Repair

- HDAC Fluorometric & Colorimetric Assays & Drug Discovery Kits
- HAT Colorimetric Assay Kit & Reagents
- DNA Damage Quantification Kit
- Glutathione & Nitric Oxide Fluorometric & Colorimetric Assay Kits

Signal Transduction

- cAMP & cGMP Assay Kits
- Akt & JNK Activity Assay Kits
- Beta-Secretase Activity Assay Kit

Adipocyte & Lipid Transfer

- Recombinant Adiponectin, Survivin, & Leptin
- CETP Activity Assay & Drug Discovery Kits
- PLTP Activity Assay & Drug Discovery Kits
- Total Cholesterol Quantification Kit

Molecular Biology & Reporter Assays

- siRNA Vectors
- Cloning Insert Quick Screening Kit
- Mitochondrial & Genomic DNA Isolation Kits
- 5 Minutes DNA Ligation Kit
- 20 Minutes Gel Staining/Destaining Kit
- β-Galactosidase Staining Kit & Luciferase Reporter Assay Kit

Growth Factors and Cytokines

Monoclonal and Polyclonal Antibodies