

**Purified Mouse Anti-human/mouse/rabbit
Actin (muscle) Antibody *HHF35,
monoclonal***

Catalog number: V103030

Unit size: 0.1 mg

Product Details

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	Lot specific (please consult certificate of analysis for given lot)
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 15 mM sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human, mouse, rabbit, chicken, non-human primates, dog, rat, cat
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Actin (muscle)
Clone	HHF35

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	WB, IHC(P), IHC(F)

Applications

Actin, aortic smooth muscle (sometimes called Cell growth-inhibiting gene 46 protein or α -actin-2) is a protein with a molecular weight of 49 kDa, found in the extracellular exosome, cytosol and protein-containing complex of cells. In Homo sapiens, actin, aortic smooth muscle binds to protein kinase and ATP. It is thought to be essential to blood pressure. Actin, aortic smooth muscle promotes DNA-templated transcription and gene expression. Actin, aortic smooth muscle plays an important role in organismal processes, in particular, glomerular mesangial cell development, vascular smooth muscle contraction and response to virus. Mutations and abnormalities in actin, aortic smooth muscle have been closely linked to a number of diseases, for instance, moyamoya disease 5 (MYMY5), aortic aneurysm, familial thoracic 6 (AAT6) and multisystemic smooth muscle dysfunction syndrome (MSMDYS). Multisystemic smooth muscle dysfunction syndrome, an autosomal dominant inheritance disorder, characterized by dilatation of the cerebral artery, volvulus of midgut and hyperperistalsis, has especially been of interest to investigators.