VHL monoclonal antibody (AT82B10)

Von Hippel-Lindau disease (VHL) is a dominant inherited syndrome characterized by the predisposition to develop various kinds of benign and malignant tumors. VHL syndrome is caused by germline mutation in the VHL tumor suppressor, and VHL tumors are associated with loss or mutation of the remaining wild-type allele. VHL protein is also involved in the degradation of hypoxia-inducible factor (HIF).

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS746-0100

100µl

Manuals, SDS & CofA

View Online »

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Von Hippel-Lindau tumor suppressor isoform 1

Application ELISA, ICC, IF, WB

Clone AT82B10

Formulation Liquid. In PBS, pH 7.4, containing 0.02% sodium azide

and 10% glycerol.

Host Mouse

Immunogen Recombinant human VHL (aa 1-154) purified from E. coli.

lgG2bk

Purity Detail Protein A affinity purified

Source Purified from ascites.

Species Reactivity Human

UniProt ID P40337

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

