

PAX6 antibody (pAb)

Catalog Nos: 61611, 61612

RRID: AB_2793701

Isotype: IgG

Application(s): WB

Reactivity: Human

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 52 kDa

Background: PAX6 (Paired Box 6) is a transcription factor with important functions in the development of the eye, nose, central nervous system and pancreas.

Required for the differentiation of pancreatic islet alpha cells (By similarity).

Competes with PAX4 in binding to a common element in the glucagon, insulin and somatostatin promoters. Regulates specification of the ventral neuron subtypes by establishing the correct progenitor domains (By similarity). Isoform 5a appears to function as a molecular switch that specifies target genes.

Immunogen: This antibody was raised against a peptide within the C-terminal region of human PAX6.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

WB*: 1:500 - 1:1,000 dilution

*Note: many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western blot.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years.

Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.

— 195

— 142

— 96

— 71

— 48

— 33

— 28

PAX6 antibody (pAb) tested by Western blot.

Detection of PAX6 by Western blot. The analysis was performed using 30 μ g HEK293 nuclear extract and PAX6 (pAb) at a 1:500 dilution.