

## PDX1 antibody (pAb)

**Catalog Nos:** 61289, 61290

**RRID:** AB\_2793583

**Isotype:** IgG

**Application(s):** WB

**Reactivity:** Human

**Volumes:** 100  $\mu$ l, 10  $\mu$ l

**Purification:** Affinity Purified

**Host:** Rabbit

**Molecular Weight:** 42 kDa

**Background:** PDX1 (pancreatic and duodenal homeobox 1, GSF, IPF1, IUF1, IDX-1, MODY4, PDX-1, STF-1) activates insulin, somatostatin, glucokinase, islet amyloid polypeptide and glucose transporter type 2 gene transcription. Particularly involved in glucose-dependent regulation of insulin gene transcription. Binds preferentially the DNA motif 5'-[CT]TAAT[ TG]-3'. During development, specifies the early pancreatic epithelium, permitting its proliferation, branching and subsequent differentiation. At adult stage, required for maintaining the hormone-producing phenotype of the beta-cell.

**Immunogen:** This PDX1 antibody was raised against a peptide within the N-terminal region of human PDX1.

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

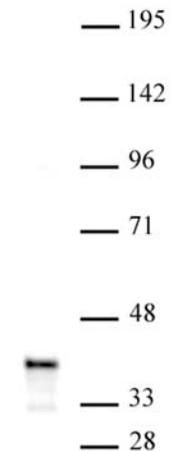
### Application Notes:

Applications Validated by Active Motif:

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

**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.



**PDX1 antibody (pAb) tested by Western blot.**  
PDX1 detection by Western blot. The analysis was performed using 30  $\mu$ g Caco-2 nuclear cell extract and PDX1 at a 1:500 dilution.