PAX7 monoclonal antibody (PBPX7)

The Pax gene family of nuclear transcription factors is comprised of nine members that function during embryogenesis to regulate the temporal and position-dependent differentiation of cells. Pax-7 is a protein specifically expressed in cultured satellite cell-derived myoblasts. In situ hybridization reveals that Pax-7 is also expressed in satellite cells residing in adult muscle. A chromosomal aberration in the gene encoding Pax-7 causes rhabdomyosarcoma 2 (RMS2).

This antibody is covered by our Worry-Free Guarantee.

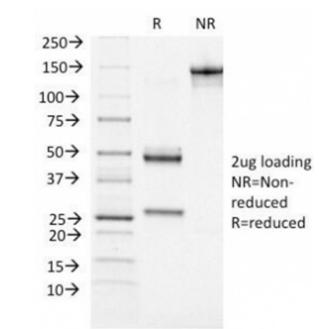
Ordering Information

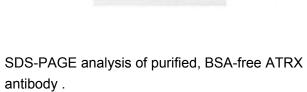
Order Online »

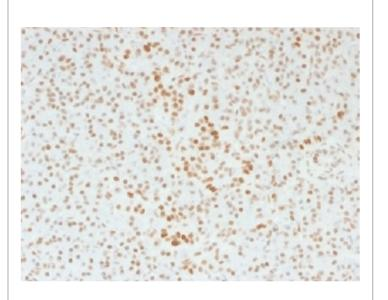
ENZ-ABS791-0100 100µg

Manuals, SDS & CofA

View Online »







IHC testing of FFPE human pancreas with ATRX antibody. Required HIER: steam section in pH6 citrate buffer for 20 min and allow to cool prior to staining.

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 Pairex box protein Pax-7, HuP1

Application IHC, WB

Clone PBPX7

Formulation Liquid. In 1X PBS containing 0.1 mg/ml BSA and 0.05%

sodium azide.

Host Mouse

Immunogen Recombinant PAX7 fragment (aa 300-600).

Isotype IgG1ĸ

Purity Detail Protein G affinity purified.

Species Reactivity Human

UniProt ID P23759

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

.

Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 info-

uk@enzolifesciences.com