# NeuroD1 polyclonal antibody

The anti-NEUROD1 antibody is designed to detect human NEUROD1 (Neuronal Differentiation 1, also known as BETA2).

Loss of NEUROD1 function has been linked to a range of disorders, including ataxia, cerebellar hypoplasia, sensorineural deafness, and severe retinal dystrophy in mice. In humans, mutations or dysregulation of NEUROD1 are associated with diseases such as Maturity-Onset Diabetes of the Young (MODY6) and noninsulin-dependent diabetes mellitus (type 2 diabetes). This dual role in both endocrine and neuronal systems makes NEUROD1 a critical marker for studying developmental biology, diabetes, and neurodegeneration.

This antibody is covered by our Worry-Free Guarantee.

# **Ordering Information**

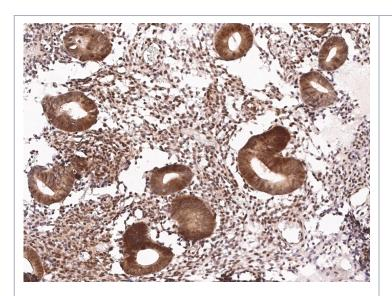
**Order Online** »

ENZ-ABS1044-0100

100µl

Manuals, SDS & CofA

**View Online** »



Immunohistochemical staining of human NEUROD1 in human endometrium with rabbit polyclonal antibody (1:2000, formalin-fixed paraffin embedded sections).

## **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 BETA2, Neurogenic differentiation factor 1

**Application** IHC (PS)

Formulation Liquid. 0.2 µm filtered solution in PBS. Carrier-free.

Preservative-free.

**Host** Rabbit

**Immunogen** Synthetic peptide corresponding to the N-terminus of

human NeuroD1.

**Isotype** IgG

Purity Detail Protein A and antigen affinity purified.

**Recommendation Dilutions/Conditions** Immunohistochemistry (paraffin sections) (1:1'000-1:4'000)

Suggested dilutions/conditions may not be available for all

applications.

Optimal conditions must be determined individually for

each application.

Species Reactivity Human

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

•



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
infouk@enzolifesciences.com