

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](#).

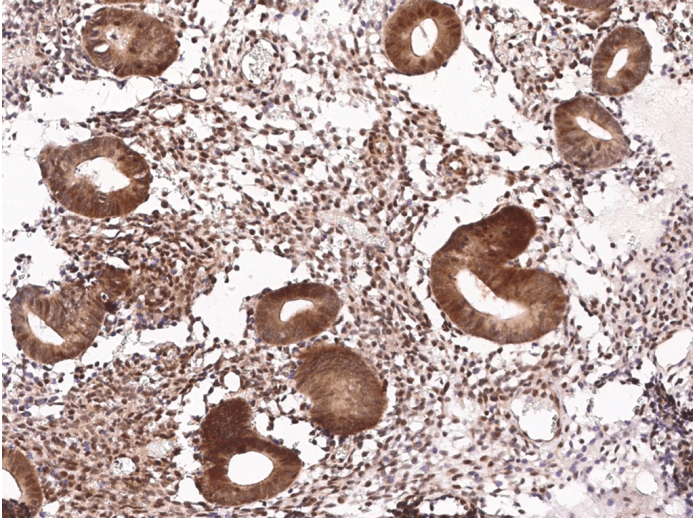
Ordering Information

[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 Immunohistochemistry (paraffin sections) (1:1'000-1:4'000)

Dilutions/Conditions 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](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com