

# NKX2.2 Recombinant monoclonal antibody (NX2/1422R)

NKX2.2 acts as a transcriptional activator. Expression of NKX2.2 has been found in neuroendocrine tumors of the gut, making it a potential marker for the study of gastrointestinal neuroendocrine tumors. More recently, NKX2.2 protein was identified as a target of EWS-FLI-1, the fusion protein specific to Ewing sarcoma, and was shown to be differentially upregulated in Ewing sarcoma on the basis of array-based gene expression analysis.

This antibody is covered by our [Worry-Free Guarantee](#).

## Ordering Information

[Order Online »](#)

ENZ-ABS370-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Homeobox protein Nkx-2.2

**Application** Flow Cytometry, IF, IHC (PS)

**Application Notes** Antibody to NKX2.2 detects a valuable marker for Ewing sarcoma, with a sensitivity of 93% and a specificity of 89%, and aids in the differential diagnosis of small round cell tumors.

**Clone** NX2/1422R

**Formulation** Liquid. In 1X PBS containing 0.1 mg/ml BSA (US sourced) and 0.05% sodium azide.

**Gene/Protein Identifier** 4821 (Gene ID)

**Host** Rabbit

**Immunogen** Recombinant human NKX2.2 protein fragment.

**Isotype** IgGk

**Purity Detail** Protein A affinity purified.

**Recommendation** Flow Cytometry (0.5-1 µg/106 cells in 0.1 mL)Immunofluorescence (1-2 µg/mL)Immunohistochemistry (formalin-fixed paraffin sections) (1-2 µg/mL for 30 minutes at RT)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

**Species Reactivity** Human

**UniProt ID** O95096



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

European Sales Office  
ENZO LIFE SCIENCES  
(ELSI) AS  
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

This antibody is covered by our **Worry-Free Guarantee**.