

# Globo H (human) monoclonal antibody (VK9)

Globo H is a hexasaccharide formed from the pentasaccharide precursor, SSEA3 (stage-specific embryonic antigen 3). It is a member of a family of antigenic carbohydrates expressed on the surface of cancer cells as glycolipids and possibly glycoproteins. Globo H is overexpressed on many carcinomas, such as breast, colon, ovarian, gastric, pancreatic, lung, and prostate. Expression of Globo H in carcinomas include more than half of all ductal, lobular, and tubular breast carcinomas, but not in non-epithelial breast tumors. Cancer stem cells and human embryonic stem cells also express global H, although expression decreases when cells are differentiated toward both neural progenitor cells and to a lesser extent in endodermal lineage cells. Expression in normal tissues is limited to cells composing secretory borders. Globo H is considered a target for immunotherapy of many epithelial cancers.

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

## Ordering Information

[Order Online »](#)

ENZ-ABS379-0050	50µg
-----------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Application** Flow Cytometry, Fluorescence microscopy, IHC (PS), WB

**Clone** VK9

**Formulation** Liquid. 0.2 µM filtered solution in PBS, pH 7.2, containing 0.09% sodium azide.

**Host** Mouse

**Immunogen** Globo H-KLH (keyhole limpet hemocyanin) synthetic conjugate

**Isotype** IgG3κ

**Purity Detail** Affinity chromatography purified.

**Recommendation** Western Blot (5 µg/mL) Immunohistochemistry (paraffin sections, 5 µg/mL) Suggested  
**Dilutions/Conditions** 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](mailto:info-usa@enzolifesciences.com)

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

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

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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