

EUROPE

KANEKA EUROGENTEC SA - LIEGE SCIENCE PARK • 4102 Seraing • BELGIUM • Tel.: +32 4 372 74 00 •
Fax: +32 4 372 75 00 Toll-free: + 800 666 00 123 • info@eurogentec.com • www.eurogentec.com

NORTH AMERICA

ANASPEC - 34801 Campus Drive • Fremont, CA 94555 • USA • Tel.: +1-510-791-9560
Toll-free: +1 800-452-5530 • Fax: +1 510-791-9572 • service@anaspec.com • www.anaspec.com

Kaneka Eurogentec products are sold for research or laboratory use only and are not to be administered to humans or used for medical diagnostics.

Technical Data Sheet

Rabbit anti-VHH, polyclonal, HiLyte™ Fluor 488 labeled EG-VHH001-HF488

100 µg
0.5 mg/mL in PBS + 0.05% NaN₃

Host: Rabbit

Isotype: IgG

Reactivity: Llama VHHs from different framework sequence families¹

Label: HiLyte™ Fluor 488

Antibody description

Besides the conventional IgGs (IgG1, composed of heavy and light chains), llamas and other camelids produce single chain IgGs (IgG2 and IgG3). They are composed of heavy chains only and are lacking the light chains and the CH1 domains. The antigen binding domain of IgG2 and IgG3, called V_HH, is the smallest naturally occurring antibody binding domain.

This secondary antibody is broad VHH-specific. It allows detecting VHH's from multiple germline sequence families. This antibody was purified using protein A.

It is conjugated to HiLyte™ Fluor 488, a green fluorophore with brighter fluorescence and better photostability than fluoresceins such as FITC. The fluorescence of HiLyte™ Fluor 488 can be observed at excitation/emission wavelength of 499 nm/523 nm. The degree of substitution (DOS) of this antibody conjugate is batch dependent and reported on the certificate of analysis.

Storage conditions

Store at 4°C.

Application

The antibody has been tested in ELISA, immunohistochemistry, immunoblotting (WB) and other immunodetection methods.

Suggested Working Dilution²:

ELISA: 5000x to 10000x

WB: 2000x

Immuno-histochemistry: 2000x

¹ We do not exclude that this secondary Ab also binds to VHH from other camelidae

² The optimal antibody dilution should be determined for each specific assay condition

For further information please contact our Customer Help Desk

For Europe:

E-mail: info@eurogentec.com

Tel: +32 4 372 76 65 • Toll-free: + 800 666 00 123

For USA:

E-mail: service@anaspec.com

Tel.: +1-510-791-9560 • Toll-free: +1-800-452-5530