

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 EG-VHH001-0025

25 µg

1 mg/mL in PBS + 0.03 % Thimerosal

Host: Rabbit

Isotype: IgG

Reactivity: Llama VHHs from different framework sequence families¹

Label: None

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 polyclonal secondary antibody generated using VHHs from multiple sources allows detection of a very large panel from different germline sequence families VHHs. This antibody was purified using protein A.

Storage conditions

Divide the solution into aliquots and store at -20°C or add an equal volume of glycerol (ACS grade or higher) and store at -20°C without aliquoting. Avoid multiple thaw-freeze cycles.

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 camelidaes

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