

Goat anti-rabbit IgG, polyclonal antibody (AP conjugate)

This antibody conjugate is used in the AMP'D[®] HSP70 high sensitivity ELISA kit (Prod. no. ENZ-KIT-101) and has been successfully employed for use with the APP Δ C31 ELISA kit (Prod. no. ADI-900-227) as part of our AMP'D[®] detection system to increase the sensitivity of this ELISA ~10-fold.

Immunoglobulin G (IgG) is an immunoglobulin monomer consisting of two (gamma) heavy chains and two light chains. Each IgG molecule contains two antigen binding domains and a single effector (Fc) domain.

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

Citations: 7

[View Online »](#)

Ordering Information

[Order Online »](#)

| | |
|-----------------|-----|
| ENZ-ABS257-5000 | 5ml |
|-----------------|-----|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ELISA

Application Notes For certified AMP'D[®] ELISA kit development.

Formulation Liquid (100 ng/ml). In 0.05M TRIS-HCl, pH 8.0, containing 0.075M sodium chloride, 0.05% hydorol M, 0.5% BSA, 0.05% Tween-20, and 50% Stabilzyme-AP.

Host Goat

Immunogen Purified rabbit IgG (whole molecule).

Purity Detail Immunoaffinity purified.

Quality Control Goat antiserum immunoaffintity purified to isolate specific anti-rabbit IgG antibodies then adsorbed with human IgG and conjugated to alkaline phosphatase.

Source Purified from rabbit serum.

Species Reactivity Rabbit

Specificity Recognizes rabbit IgG.

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)