

Head-Peptide polyclonal antibody

Ordering Information

[Order Online »](#)

| | |
|-------------|------|
| JBS-ABD-058 | 20µl |
|-------------|------|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

| | |
|-------------------|---------------------------|
| Handling | Avoid freeze/thaw cycles. |
| Long Term Storage | -20°C |
| Shipping | Blue Ice |

Regulatory Status RUO - Research Use Only

Product Details

| | |
|--------------------------------|---|
| Application | ELISA, ICC |
| Formulation | liquid (Supplied as serum, preserved in glycerol) |
| Immunogen | Raised against the Periplaneta americana midgut neuropeptide, ANRSPSLRLRLRFamide, coupled to bovine thyroglobulin using glutaraldehyde. |
| Technical Info / Product Notes | For the Original Manufacturer's data sheet please click here . |



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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
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