

# IL-34 Antibody

## Ordering Information

[Order Online »](#)

|                |       |
|----------------|-------|
| FGX-IL34-101AP | 100µg |
|----------------|-------|

## Manuals, SDS & CofA

[View Online »](#)

# Handling & Storage

**Regulatory Status** RUO - Research Use Only

## Product Details

|                           |                   |
|---------------------------|-------------------|
| <b>Application</b>        | ELISA, IP, WB     |
| <b>Host</b>               | Rabbit            |
| <b>Immunogen</b>          | Interleukin-34    |
| <b>Isotype</b>            | IgG               |
| <b>Purity Detail</b>      | Affinity Purified |
| <b>Species Reactivity</b> | Human, Mouse, Rat |
| <b>UniProt ID</b>         | Q6ZMJ4            |



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)