

Anti-Human Interferon Lambda Receptor 1, Clone HLR-14 (MAb)

Ordering Information

[Order Online »](#)

| | |
|-------------|------|
| PBL-21886-1 | 50µg |
|-------------|------|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

| | |
|---------------------------|---------------------------|
| Handling | Avoid freeze/thaw cycles. |
| Short Term Storage | -20°C |
| Long Term Storage | -80°C |
| Shipping | Dry Ice |

Regulatory Status RUO - Research Use Only

Product Details

| | |
|---------------------------|---|
| Alternative Name | IFN-λ, Interferon-λ |
| Application | ELISA, Flow Cytometry |
| Clone | HLR14 |
| Formulation | 0.3 M Sodium Bicarbonate, 0.2 M Sodium Chloride, pH 8.0 |
| Host | Mouse |
| Isotype | IgG1 |
| Purity | > 95% |
| Purity Detail | Protein G purified. |
| Species Reactivity | Human |
| Specificity | Binds to extracellular domains of human interferon lambda receptor. |



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)