

Human IL-1 β ELISA Standard Recombinant Protein, eBioscience™

Catalog Number 29-8018-65

Pub. No. MAN0018027 Rev. A.0 (30)

 **WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from thermofisher.com/support.

Product description

| | |
|---------------------------|--|
| Contents | Human IL-1 β ELISA Standard Recombinant Protein, eBioscience™ |
| Size | 4 pack |
| Source | <i>E. coli</i> expressed amino acids Ala117-Ser269 of human IL-1 beta (accession # NM_000576). |
| Formulation | Lyophilized from aqueous buffer, may contain carrier protein/stabilizer |
| Handling | <p>Reconstitute lyophilized standard using distilled water. For reconstitution volume see vial label. Allow the standard to reconstitute for 10-30 minutes. Swirl or mix gently to ensure complete and homogeneous solubilization. Concentration of reconstituted standard is 150 pg/mL.</p> <p>Note: Single-use vials. Use within 4 hours of reconstitution and discard after use. Dilute only as shown above to ensure accurate recovery.</p> |
| Storage Conditions | Store at 2-8°C. |

thermofisher.com/support | thermofisher.com/askaquestion



Manufacturer: Bender MedSystems GmbH | Campus Vienna Biocenter 2 | 1030 Vienna, Austria

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.