


Rat IgG2a ELISA Standard Protein, eBioscience™

Catalog Number 39-50510-65

Pub. No. MAN0018155 Rev. B.0 (31)

 **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](https://www.thermofisher.com/support).

Product description

Contents	Rat IgG2a ELISA Standard Protein , eBioscience™
Size	4 pack
Formulation	Lyophilized from aqueous buffer, may contain carrier protein/stabilizer
Handling	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 50 ng/mL. Note: Mix well prior to making dilutions. The standard vial has to be used immediately after reconstitution and cannot be stored.
Storage Conditions	Store at 2-8°C.

[thermofisher.com/support](https://www.thermofisher.com/support) | [thermofisher.com/askaquestion](https://www.thermofisher.com/askaquestion)



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

For descriptions of symbols on product labels or product documents, go to [thermofisher.com/symbols-definition](https://www.thermofisher.com/symbols-definition).

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.

©2021 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.