



MONOCLONAL ANTIBODY ANTI-IgE

Product Data Sheet

Part # MABETS07 Version : 2

For research use only. Not for use in diagnostic procedures.

PRODUCT SPECIFICATIONS

Description: Mouse monoclonal ETS07 anti-human IgE antibody

Host species: Mouse

Form: Lyophilized

Purity: Protein A affinity purified

Buffer: P04 50 mM pH7, 1% BSA, 1% trehalose, 0.5% mannitol

Storage: 2° - 8°C.

Once reconstituted, the antibody solution may be frozen at -60°C or below, and can be thawed only once. To avoid freeze/thaw cycles it is recommended to dispense remaining antibody solution into disposable plastic vials for storage at -60°C or below

Specificity: This antibody reacts with heavy chains on IgE but not with other immunoglobulins in human serum.

RECONSTITUTION

Reconstitute with 0.5 mL of sterile distilled water. The antibody concentration will be 0.2 mg/mL.

The reconstitution volume can be adjusted from 0.3 mL to 1 mL depending on the application.

APPLICATIONS

- The monoclonal antibody anti-IgE has been validated in a functional assay to monitor the histamine release upon treatment of fresh blood donors. Please refer to the application note "Rapid evaluation of compound allergenicity by measuring histamine release on stimulated whole blood from different donors" for in-depth information.
- The monoclonal antibody anti-IgE can be used for the detection of total IgE level and IgE antibodies.



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF

Manufactured by Cisbio Bioassays - Parc Marcel Boiteux - 30200 Codolet - FRANCE

www.revvity.com

revvity

Revvity, Inc.
940 Winter Street
Waltham, MA 02451 USA
www.revvity.com

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2023, Revvity, Inc. All rights reserved.