



G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

WellCoat™ Stabilizer

(Stabilizer for ELISA Coating Plate)

(Cat. # 786-1217)



think proteins! think G-Biosciences www.GBiosciences.com

INTRODUCTION

WellCoat™ Stabilizer is developed to be used as a stabilizing agent for antigens, antibodies or other ligands coated on microwell ELISA plates. WellCoat™ Stabilizer created a protective layer over the immobilized agents, allowing storage of the coated plates for an extended period of time without significant loss of signal. The stabilizer may be used directly to coat the pretreated plates or mix with a blocking agent- suitable for any buffer combination. Use of WellCoat™ Stabilizer eliminates washing steps prior to storage of the coated plates - offer considerable time saving when preparing a large number for pre-coated plates for storage.

ITEM(S) SUPPLIED

Cat. #	Description	Size
786-1217	WellCoat™ Stabilizer	1L

STORAGE CONDITIONS

WellCoat™ Stabilizer is shipped at ambient temperature. Upon arrival, store it at 4°C. If stored properly as instructed, the product is stable for 1 Year.

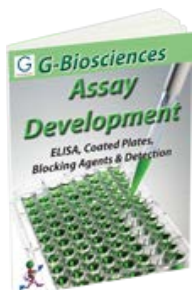
ADDITIONAL ITEMS REQUIRED

- Blocking agent.
- ELISA plate sealing covers or pouches.
- Desiccant pouches

PROTOCOL

1. After the treatment with blocking agent aspirate and remove the blocking solution.
2. Apply 300 µl (fill the wells) with WellCoat™ Stabilizer. Alternatively, mix WellCoat™ Stabilizer with blocking agent in 1:1 ratio and then apply to the ELISA microwells.
3. Let the ELISA plate stand for 30 minutes at Room Temperature.
4. Gently aspirate off the contents of the ELISA microwells and allow the plate to air dry at Room Temperature.
5. Seal the ELISA plate with a pouch containing desiccant.
6. Store the plates at 4°C for future use.
NOTE: *The Wellcoat™ Stabilizer treated ELISA plates/microwells can be stored at 4°C for 1 year without significant loss of signal.*
7. *Before use, wash the microwells three times with PBS*

RELATED PRODUCTS



Download our Assay Development Handbook <http://info2.gbiosciences.com/complete-assay-development-handbook>

For other related products, visit our website at www.GBiosciences.com or contact us.

Last saved: 1/25/2017 CMH



www.GBiosciences.com