



052PR-02

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

A Geno Technology, Inc. (USA) brand name

Well-Coated™ Glutathione

96-Well Plates Coated with Glutathione
for Binding GST Tagged Proteins

(Cat. # 786-751, 786-764, 786-765)



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

INTRODUCTION

Well-Coated™ Glutathione plates are designed to specifically bind GST (Glutathione S-Transferase) tagged proteins and peptides. The plates have immobilized glutathione and can be used for the isolation of GST-tagged proteins direct from bacterial lysates and then used for subsequent ELISA protocols.

The wells are coated to a 100µl depth and are supplied pre-blocked in our proprietary Superior™ Blocking Buffer. The plates are protected with our WellCoat™ Stabilizer (Cat. # 786-1217) that creates a protective layer over the immobilized agents. The reagent will not interfere with the assay and has no effect of the efficiency or capacity of the wells. The WellCoat™ Stabilizer offers greater protection and shelf life of the plates. In some cases, the protective layer may give the appearance of a white coating. The clear, white and black plates are offered for colorimetric, chemiluminescence and fluorescent detection systems, respectively.

ITEM(S) SUPPLIED

Cat. #	Components	Size
786-751	Well-Coated™ Glutathione 8-well strip plate, Clear	5 plates
786-764	Well-Coated™ Glutathione 96 well plate, Black	5 plates
786-765	Well-Coated™ Glutathione 96 well plate, White	5 plates

STORAGE CONDITIONS

Shipped at ambient temperature. Upon arrival, store unopened at 4°C. Once opened the plates can be stored in a resealable bag (ZipLoc) with an appropriate desiccant at 4°C.

BINDING CAPACITY

Well-Coated™ Glutathione: ~10ng purified GST/well

PROTOCOL

The following protocol is a simple direct ELISA protocol and the protocol and reagents used will have to be optimized for specific applications and assays.

Additional Item(s) Required

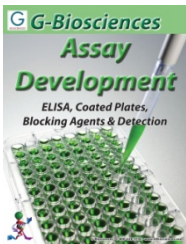
- GST tagged protein cell lysate
- Dilution Buffer: Tris buffered saline (TBS, Cat. # 786-288) or phosphate buffered saline (PBS, Cat. # 786-289).
- Wash Buffer: femtoTBST™ (Cat. # 786-161) or femtoPBST™ (Cat. # 786-162); 10X concentrated wash buffers supplemented with Tween® 20. Alternatively use PBS supplemented with 0.05% Tween® 20.
- Enzyme Labeled Primary Antibody against the GST tagged protein; visit www.GBiosciences.com for horseradish peroxidase (HRP) and alkaline phosphatase (AP) labeling kits.
- Detection system, femtoELISA™ is a chromogenic detection system for HRP and AP (Cat. # 786-110 to 786-113)

Direct ELISA Assay

1. Wash the wells to be used three times with 200µl Wash Buffer.
2. Dilute the cell lysate with Dilution Buffer and add up to 100µl diluted lysate to each well.
3. Incubate at room temperature for 1-2 hours, for optimal binding use a plate shaker.
4. Wash each well three times with 200µl Wash Buffer.
5. Add 100µl enzyme labeled primary antibody.
6. Incubate at room temperature for 0.5-1 hour with shaking.
7. Wash each well three times with 200µl Wash Buffer.
8. Detect the label signal according to the manufacturer's instructions, using 100µl detection reagent per well.

RELATED PRODUCTS

Download our Assay Development Handbook.



<http://info.gbiosciences.com/complete-assay-development-handbook>

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



www.GBiosciences.com