

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

ELISA Plate Coating Buffer (5X)

Catalog #: 4500-050-01

Size: 50 ml

Catalog #: 4500-100-01

Size: 100 ml

Catalog #: 4500-01k-01

Size: 1 liter

Background: Trevigen ELISA Plate Coating Buffer, with an optimal formulation of buffers and salts, can improve assay performance by maximizing the ELISA signal and decreasing the background noise, or increase the signal to noise ratio. Using Trevigen ELISA Plate Coating Buffer can also extend shelf life time of the coated plates.

Applications: Dilute 5X Coating Buffer to 1X with deionized water prior for use. Simply add your protein, antigen or antibody to the diluted 1X working coating buffer, mix and pipette onto the plate.

Storage: 4 °C storage.

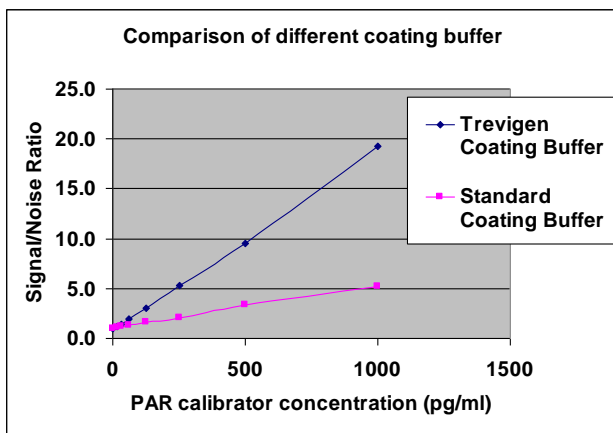


Fig. 1 Comparison of ELISA performance by different coating buffer.

In one example, Trevigen ELISA plate coating buffer increases the ELISA Performance by increasing the signal to noise ratio 4-5 times more than normal coating buffer. Standard coating buffer: Sigma-Aldrich Carbonate-bicarbonate buffer (Cat# C3041-50CAP).

Related Products:

Catalog #	Description	Size
3455-096-03	10X Coating Buffer	1 ml
4671-096-02	20X PARP Buffer	500 µl
4830-250-02	10X Binding Buffer	20 ml

ELISA Plate Coating Buffer

Catalog #: 4500-050-01,
4500-100-01, 4500-01k-01

Storage: 4 °C

TREVIGEN®
1-800-873-8443