

Solid Phase Blocking Buffer, 10X

Catalog Number X109-25ML

Catalog Number X109-250ML



ARBOR
ASSAYS

FEATURES

- Blocking buffer for solid phases
- 10X Concentrate
- Stable solution at 4°C

INTRODUCTION

Solid phase immunoassays utilizing microtiter plates are the predominant assay format for the detection of low concentration biomolecules in biological samples. Typical formats include EIA, ELISA, CIA and CLIA type assays where either an antigen or an antibody is immobilized onto a high binding solid phase surface. Detection of the specific biomolecule utilize either labeled antigen or antibody and utilize enzymes such as HRP to give rise to a colorimetric or chemiluminescent signal.

Both antigen and antibody coated plates can be generated using a suitable coating buffer and after being blocked and dried these plates are stable for many years. The Arbor Assays Solid Phase Blocking Buffer is optimized to block uncoated sites after the attachment of antigens and antibodies to microtiter plates. Purified antibodies and antigens are attached to the wells of a suitable microtiter plate in a buffer such as our Coating Buffer, X108-10ML or X108-100ML. After incubation, typically overnight, the contents of the wells are discarded and the wells filled with diluted Blocking Buffer, X109-25ML or X109-250ML. After incubation, again typically overnight, the plates are dried and can be stored at 4°C for extended periods of time.

FORM:	10X Concentrate
STORAGE:	4°C
STABILITY:	Until expiration date on container
USES:	As a microtiter plate blocking buffer after attachment of antigens and antibodies
SUGGESTED USE:	Dilute one part with 9 parts of deionized water
COUNTRY OF ORIGIN:	USA

FOR RESEARCH USE ONLY