

Solid Phase Coating Buffer, 20X

Catalog Number X108-10ML

Catalog Number X108-100ML



ARBOR
ASSAYS

FEATURES

- Coating buffer for solid phase attachment of antibodies
- 20X 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 Coating Buffer is optimized to allow the attachment of antigens and antibodies to microtiter plates by passive absorption. Purified antibodies and antigens are diluted into the diluted coating buffer and added to the wells of a suitable microtiter plate. After incubation, typically overnight, the contents of the wells are discarded and the wells filled with a blocking buffer, such as our 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:	20X Concentrate
STORAGE:	4°C
STABILITY:	Until expiration date on container
USES:	As a microtiter plate coating buffer for attachment of antigens and antibodies
SUGGESTED USE:	Dilute one part with 19 parts of deionized water
COUNTRY OF ORIGIN:	USA

FOR RESEARCH USE ONLY