

MESO SCALE DISCOVERY® Assay Reagents

MSD® Blocker A Kits

Buffer	Part Number	Quantity
MSD Blocker A Kit, 1L	R93AA-1	1L
MSD Phosphate Buffer (5X)	R93SA-1*	200 mL
MSD Blocker A	R93BA-1*	50 g
MSD Blocker A Kit, 250 mL	R93AA-2	250 mL
MSD Phosphate Buffer (5X)	R93SA-2*	50 mL
MSD Blocker A	R93BA-2*	12.5 g

* These items are not sold separately

Storage	Store MSD Phosphate Buffer (5X) at room temperature. Store reconstituted MSD Blocker A Solution at 4°C.
Stability	When stored properly, MSD Phosphate Buffer (5X) and MSD Blocker A (dry powder) are stable until the expiration date printed on the product label. Once prepared, MSD Blocker A Solution is stable for 5 weeks after the date of reconstitution.
Description	MSD Blocker A Solution is a cocktail of proteins in a PBS-based buffer optimized for use with MSD MULTI-ARRAY® and MULTI-SPOT® Assays. MSD Blocker A Solution "blocks" non-specific binding of proteins to the plate surface to enhance sensitivity and reduce background in some assays.
Preparation Procedure for MSD Blocker A Solution	<p>Step 1: Addition of deionized water to the MSD Blocker A dry powder. Add 800 mL of deionized water to the 1 L bottle. OR Add 200 mL of deionized water to the 250 mL bottle.</p> <p>Step 2: Soak and stir. Allow to soak for 2 hours. Add a magnetic stir bar and stir until all protein is resuspended (30 min. - 2 hrs.). Upon dissolution of all material, the solution will be pale yellow and clear.</p> <p>Step 3: Add MSD Phosphate Buffer (5X). Add entire contents of MSD Phosphate Buffer (5X) bottle to the reconstituted solution (200 mL for the 1 L bottle, 50 mL for the 250 mL bottle and stir for an additional 10 minutes. Vacuum filter the prepared MSD Blocker A Solution with a 0.2 µm filter into a clean storage container. Attach the separate MSD Blocker A Solution label (provided) to the new container. Store at 4°C. <i>Following the instructions listed above, the MSD Blocker A Solution is a 5% solution of protein in a PBS-based buffer.</i></p>
Additional Notes	<p>To prepare Blocker A Solution for one plate use only, follow steps indicated below: Weigh out 1.25 g of Blocker A dry powder Add 20 mL of deionized water Stir until all protein is resuspended Add 5 mL of MSD Phosphate Buffer (5X)</p> <p>A 5% MSD Blocker A Solution may be prepared in buffers other than MSD Phosphate Buffer (5X), depending upon the research application. For example, when assaying phosphoproteins, the use of a Tris-based buffer system is recommended. Follow the instructions listed above if using a 5X stock of buffer, or simply add a 1X solution of the desired buffer (250 mL or 1L as appropriate) to the Blocker A dry powder in place of a 5X buffer stock and deionized water.</p>

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC OR THERAPEUTIC PROCEDURES.

