

Meso Scale Discovery® Phosphoprotein Reagent Kit



Reagent	Quantity	Storage
Tris Wash Buffer (10X)	1L	2-8°C
Tris Lysis Buffer (1X)	200 mL	2-8°C
Phosphatase Inhibitor I (100X)	2 mL	Aliquot (100 µL) and store at 2-8°C
Phosphatase Inhibitor II (100X)	2 mL	Aliquot (100 µL) and store at 2-8°C
Protease Inhibitor Solution (50X)	4 mL	Aliquot (200 µL) and store at -20°C
Blocker A	15 g, dry powder	RT
Blocker B	2 g, dry powder	RT
Blocker D-M	3.6 mL	-20°C
Blocker D-R	1 mL	-20°C
Blocker D-G	1 mL	-20°C
Blocker D-B	3 mL	-20°C

Reagents supplied in 20-plate quantities.

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