

Product Information

Description:	This buffer is intended for use with mammalian cells during magnetic cell isolation or single cell processing. This buffer is optimized and formulated to reduce nonspecific binding and cell clumping.
Instructions:	Dilute 5-fold with sterile distilled water. For sterile conditions, filter the buffer after dilution and use under laminar flow hood.
Quantity:	25 ml
Quality:	Prepared under sterile conditions and filtered (0.2 μ m)
Storage:	Upon receipt, the buffer should be stored at 4°C
Stability:	When stored at 4°C, the product is stable for one year