

Product Information

Description:	FcRn Recycling Dilution Buffer (pH 6) is an acidified buffer optimized for use FcRn: IgG Recycling HMEC-1 Cell Pool (BPS Bioscience #82163).
Supplied As:	Aqueous solution.
Applications:	Use in cellular assays involving IgG recycling via FcRn as dilution buffer to allow a more efficient FcRn binding to the IgGs.
Quantity:	15 ml
Storage/Stability:	Upon receipt, store at Room Temperature. Stable for 6 months from date of receipt, when stored as directed.
Instructions for Use:	This is a ready to use buffer.
Optimized for Use With:	FcRn: IgG Recycling HMEC-1 Cell Pool (BPS Bioscience #82163)

Version 040624