

## Product Information

---

<b>Description:</b>	FcRn Recycling Wash Buffer (pH 7.0-7.4) is a neutral pH buffer optimized for use FcRn: IgG Recycling HMEC-1 Cell Pool (BPS Bioscience #82163).
<b>Supplied As:</b>	Aqueous solution.
<b>Applications:</b>	Use in cellular assays involving IgG recycling via FcRn as wash buffer to remove unbound IgG molecules.
<b>Quantity:</b>	75 ml
<b>Storage/Stability:</b>	Upon receipt, store at -80°C. Stable for 6 months from date of receipt, when stored as directed.
<b>Instructions for Use:</b>	This is a ready to use buffer.
<b>Optimized for Use With:</b>	FcRn: IgG Recycling HMEC-1 Cell Pool (BPS Bioscience #82163)

*Version 042624*