

## Product Information

<b>Description:</b>	U2 Assay Buffer is an assay buffer optimized for use with the BPS Bioscience Intrachain TR-FRET Assay Kits and Ubi-Mix™ (BPS Bioscience #82185).
<b>Supplied As:</b>	Aqueous solution.
<b>Applications:</b>	Screen molecules that inhibit ubiquitin ligase activity in drug discovery high throughput screening (HTS) applications, determine compound IC <sub>50</sub> and perform real-time kinetics analyses in Intrachain TR-FRET assays.
<b>Quantity:</b>	2 x 10 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 assay buffer.
<b>Optimized for Use With:</b>	CBL-B-Driven Tyro3 Ubiquitination Intrachain TR-FRET Assay Kit (BPS Bioscience #78338) CBL-B-driven SRC ubiquitination intrachain TR-FRET Assay Kit (BPS Bioscience #78820) CBL-B-driven AXL ubiquitination intrachain TR-FRET Assay Kit (BPS Bioscience #78821) (BPS Bioscience #78823) C-CBL-driven AXL Ubiquitination Intrachain TR-FRET Assay Kit DCAF15 Intrachain TR-FRET Assay Kit (BPS Bioscience #78543) WWP1 Intrachain TR-FRET Assay Kit (BPS Bioscience #78554) MURF3 Intrachain TR-FRET Assay Kit (BPS Bioscience #78555) ChoosE3-Freedom Intrachain TR-FRET Assay Kit (BPS Bioscience #78560) ChoosE2-Freedom Intrachain TR-FRET Assay Kit (BPS Bioscience #78561) TRAF6 Intrachain TR-FRET Assay Kit (BPS Bioscience #78598) and other Intrachain TR-FRET Assay Kits (please refer to the assay datasheet to determine applicability of U2 Assay Buffer) Ubi-Mix™ (BPS Bioscience #82185)