

Product Information

Description:	This 2X KRAS Nucleotide Exchange Buffer is optimized for use with the KRAS Nucleotide Exchange Assay Kits.
Applications:	Studying enzyme kinetics and KRAS-GDP or -GTP binding status for drug discovery and high throughput targeting of KRAS.
Quantity:	10 ml
Storage:	-20°C
Stability:	Stable for at least 6 months from date of receipt when stored as directed.
Instructions for Use:	For a detailed protocol on the use of this product, please see the assay kit protocols.
Optimized for Use With:	KRAS(G12C) Nucleotide Exchange Assay Kit (BPS Bioscience, #79859) KRAS(G12D) Nucleotide Exchange Assay Kit (BPS Bioscience, #78355)