

Product Information

Description:	This 2X KRAS(G12C) Buffer is optimized for use with the KRAS(G12C) Nucleotide Exchange Assay Kit (BPS Bioscience #79859).
Applications:	Studying enzyme kinetics and KRAS(G12C)-GDP or -GTP binding status for drug discovery and HTS targeting of KRAS(G12C).
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 assay kit protocol.
Optimized for Use With:	KRAS(G12C) Nucleotide Exchange Assay Kit (BPS Bioscience #79859).