

## Product Information

<b>Description:</b>	Guanosine 5'-triphosphate sodium salt hydrate (GTP) functions as a carrier of phosphates and pyrophosphates involved in channeling chemical energy into specific biosynthetic pathways.
<b>Purity:</b>	≥95% by HPLC
<b>Applications:</b>	Useful for studying enzyme kinetics and KRAS(G12C)-GDP or -GTP binding status for drug discovery and HTS targeting of KRAS(G12C). Suitable for use with the KRAS(G12C) Nucleotide Exchange Assay Kit (BPS Bioscience #79859) or KRAS(G12C) Coupled Nucleotide Exchange Assay Kit (SOS1 mediated) (BPS Bioscience #78004).
<b>MW:</b>	523.18 Da
<b>CAS Number:</b>	36051-31-7
<b>Empirical Formula:</b>	C <sub>10</sub> H <sub>16</sub> N <sub>5</sub> O <sub>14</sub> P <sub>3</sub> · Na
<b>Supplied As:</b>	10 mM aqueous solution
<b>Quantity:</b>	100 µl (79861-1) or 500 µl (79861-2)
<b>Storage:</b>	-20°C
<b>Stability:</b>	Stable for at least 12 months from date of receipt when stored as directed. Avoid freeze/thaw cycles.
<b>Instructions for Use:</b>	For a detailed protocol on the use of this product, please see assay kit protocol.
<b>Optimized for Use With:</b>	<ul style="list-style-type: none"><li>• KRAS(G12C) Nucleotide Exchange Assay Kit (BPS Bioscience #79859).</li><li>• KRAS(G12C) Coupled Nucleotide Exchange Assay Kit (SOS1 mediated) (BPS Bioscience #78004).</li></ul>
<b>Structure:</b>	

