

Product Information

---

<b>Description:</b>	5x Raf Substrate is a substrate designed for use in B- and C-Raf (rapidly accelerated fibrosarcoma) assays.
<b>Supplied as:</b>	Aqueous solution.
<b>Formulation:</b>	TBS (Tris Buffered Saline) and 3 mM DTT
<b>Storage:</b>	Upon receipt, store at -80°C.
<b>Stability:</b>	When stored at -80°C, the product is stable for six months after first thaw. Avoid repeated freeze/thaw cycles.
<b>Application:</b>	<p>It is suitable for use in b-Raf and c-Raf assays using purified proteins. This mixture has been optimized for use:</p> <ul style="list-style-type: none"><li>• BRAF (WT) Kinase Assay Kit (BPS Bioscience #78316)</li><li>• BRAF (V600E) Kinase Assay Kit (BPS Bioscience #48688)</li><li>• c-Raf Kinase Assay Kit (BPS Bioscience #79570)</li></ul> <p>Use in other assays may require optimization.</p>