

Product Information

---

<b>Description:</b>	FAM-cAMP is a 1 mM solution of the fluorogenic PDE Substrate 1 {2'-(6-[Fluoresceinyl]aminohexylcarbamoyl)adenosine-3',5'-cyclic monophosphate}.
<b>Supplied As:</b>	Aqueous solution.
<b>Purity:</b>	>95% by HPLC.
<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>	It is suitable for use in phosphodiesterase assays using purified proteins. It has been optimized for use with BPS Bioscience assay kits for PDE1, PDE2, PDE3, PDE4, PDE7, PDE8, PDE10 and PDE11. Use in other assays may require optimization.