

## Product Information

---

<b>Description:</b>	Histone H4 peptide, amino acids 1 to 21, acetylated at the N-terminus and biotinylated on the side chain of Lys. A GG spacer has been added on the C-terminus of Lys. Optimized for use with the PRMT5 Homogeneous Assay Kit (BPS Bioscience #52052).
<b>Supplied as:</b>	400 $\mu$ M in water (200 $\mu$ l)
<b>Size:</b>	2.5 nmole (52018-1) or 80 nmole (52018-2)
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Purity:</b>	$\geq$ 90% by HPLC
<b>MW:</b>	2602.1 Da
<b>Applications:</b>	Useful as a substrate for methyltransferases or acetyltransferases and for screening small molecular inhibitors of Histone 4 (R3) methyltransferases for drug discovery.