

Product Information

Description	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).
Supplied as:	2.5 nmole: 400 μ M in water (6.5 μ l); 80 nmole: 400 μ M in water (200 μ l)
Size:	2.5 nmole (52018-1) or 80 nmole (52018-2)
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Purity:	\geq 90% by HPLC
MW:	2602.1 Da
Application:	Useful as a substrate for methyltransferases or acetyltransferases and for screening small molecular inhibitors of Histone 4 (R3) methyltransferases for drug discovery.