

co-(5'-deoxyadenosin-5'-yl)-cobinamide, f-(dihydrogen phosphate), inner salt, 3'-ester with (5,6-dimethyl-1- α -D-ribofuranosyl-1H-benzimidazole- κ N3), hydrate

Purity Specification:	≥95%			
Molecular formula:	C72H100CoN18O17P • XH2O			
CAS number:	Formula weight:	1579.58	Expiration date:	11NOV2028

Tests	Results
Mass Spec	[M+2H] ²⁺ : 790.4
TLC	Purity: 100 %
UV	λ max: 518.00 nm λ max: 263.00 nm

Gen Labucque

WWW.CAYMANCHEM.COM