



Product Code	Product Description	G-Protein Coupled	Sizes	List Price (USD)	Vector	Availability
#Z0100N	D2 Dopamine Receptor	Gi	5mL	\$295	BacMam	immediate
#Z0101N	Human M2 Muscarinic Receptor	Gi	5mL	\$295	BacMam	immediate
#Z0102N	A1 Adenosine Receptor	Gi	5mL	\$295	BacMam	immediate
#Z0103N	Succinate Receptor	Gi	5mL	\$295	BacMam	immediate
#Z0200N	$\beta$ 2 Adrenergic Receptor	Gs	5mL	\$295	BacMam	immediate
#Z0300N	Human M1 Muscarinic Receptor	Gq	5mL	\$295	BacMam	immediate
#Z0400N	D1 Dopamine Receptor	Gs/a	5mL	\$295	BacMam	immediate
#Z0401N	Calcitonin Receptor	Gs/q	5mL	\$295	BacMam	immediate