



JBS Rainbow

Set of JBS Deep Purple, JBS True Blue, JBS Xtal Green and JBS Bright Red

Cat. No.	Amount
CO-305	4 x 300 µl

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: 12 months