



Promega Points™

Frequent Buyer Program

***Promega Points Program ends September 30, 2012.**

Point Redemption Card

Place Promega Point Here	Place Promega Point Here	Place Promega Point Here
Place Promega Point Here	Place Promega Point Here	Place Promega Point Here
Place Promega Point Here	Place Promega Point Here	Place Promega Point Here
Place Promega Point Here	Place Promega Point Here	Place Promega Point Here
Place Promega Point Here	Place Promega Point Here	Place Promega Point Here
Place Promega Point Here	Place Promega Point Here	Place Promega Point Here
Place Promega Point Here	Place Promega Point Here	Place Promega Point Here
Place Promega Point Here	Place Promega Point Here	Place Promega Point Here
Place Promega Point Here	Place Promega Point Here	Place Promega Point Here
Place Promega Point Here	Place Promega Point Here	Place Promega Point Here
Place Promega Point Here	Place Promega Point Here	Place Promega Point Here
Place Promega Point Here	Place Promega Point Here	Place Promega Point Here
Place Promega Point Here	Place Promega Point Here	Place Promega Point Here
Place Promega Point Here	Place Promega Point Here	Place Promega Point Here
Place Promega Point Here	Place Promega Point Here	Place Promega Point Here
Place Promega Point Here	Place Promega Point Here	Place Promega Point Here
Place Promega Point Here	Place Promega Point Here	Place Promega Point Here

White labels = 1 point
Yellow labels = 5 points

Please allow 4 to 6 weeks for delivery.