



Gen Labucque

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM