

Item Number	Custom Genes
SQ-GENES-XXXX	Special Quote - Custom Genes