

Cell-Free Yield Comparison

Maximum yields measured for Cell Free Expression Systems

