

Reporter Dyes for the Luminex Analyzer

