

555002

Poly[9,9-didodecylfluorenyl-2,7-yleneethynylene]

