

688096
Tris[2-phenylpyridinato-C²,N]iridium(III)

