

666025
13,6-N-Sulfinylacetamidopentacene, 97%

