



Color Table Substance

---



YDL037c



YBL071C