

Catalog # 120 Plasmid pART7 PtGFP

For cloning vector pART7 and binary vector pART27 see:
Gleave AP 1992 A versatile binary vector system with a T-DNA organisational structure conducive to efficient integration of cloned DNA into the plant genome. *Plant Molecular Biology* 20: 1203-1207.

For expression of PtGFP in plants see:
Schulte A; Lorenzen I; Böttcher M; Plieth C 2006 A novel pH indicator for expression in plants *Plant Methods* 2:7 doi:10.1186/1746-4811-2-7

