

PUC18-Gluc with SS

