

PUC18-Gluc without SS

