

Suzanne M. Hibbs, Melissa Jahangiri, Gregory Wemhoff*, and Gary Davis

72478
1030