



Spring Art Showcase March 12 - June 7, 2013

AN INTERNATIONAL CELEBRATION OF

Women *through* **art**



AN INTERNATIONAL CELEBRATION OF

Women *through* art

Inspired by International Women's Day, the spring showcase features the struggles and empowerment of women. Exhibiting are **Dr. Lilian Nabulime**, a sculptor from Uganda whose work promotes awareness about HIV/AIDS; **Michael Kienitz**, a photographer whose images document female microfinance borrowers and coffee farmers in Latin America; and **Jane Madrigal**, a Chicana artist whose woodcut prints focus on revolutionary women.

March 12 - June 7, 2013

Exhibit open to the public:

M-F 8:00 a.m. - 4:00 p.m.

Groups larger than ten

by appointment only:

(608) 298-4865

Opening Reception

Tuesday, March 12, 2013 4:30-6:30pm

BioPharmaceutical Technology Center

5445 East Cheryl Parkway, Fitchburg, WI 53711



2800 Woods Hollow Rd.
Madison, WI 53711

This exhibit is made possible through support from Promega Corporation. For future art show announcements, submit your name and email address to: artshow@promega.com

Show produced by Daniel Swadener: swadener@gmail.com | danielswadener.com