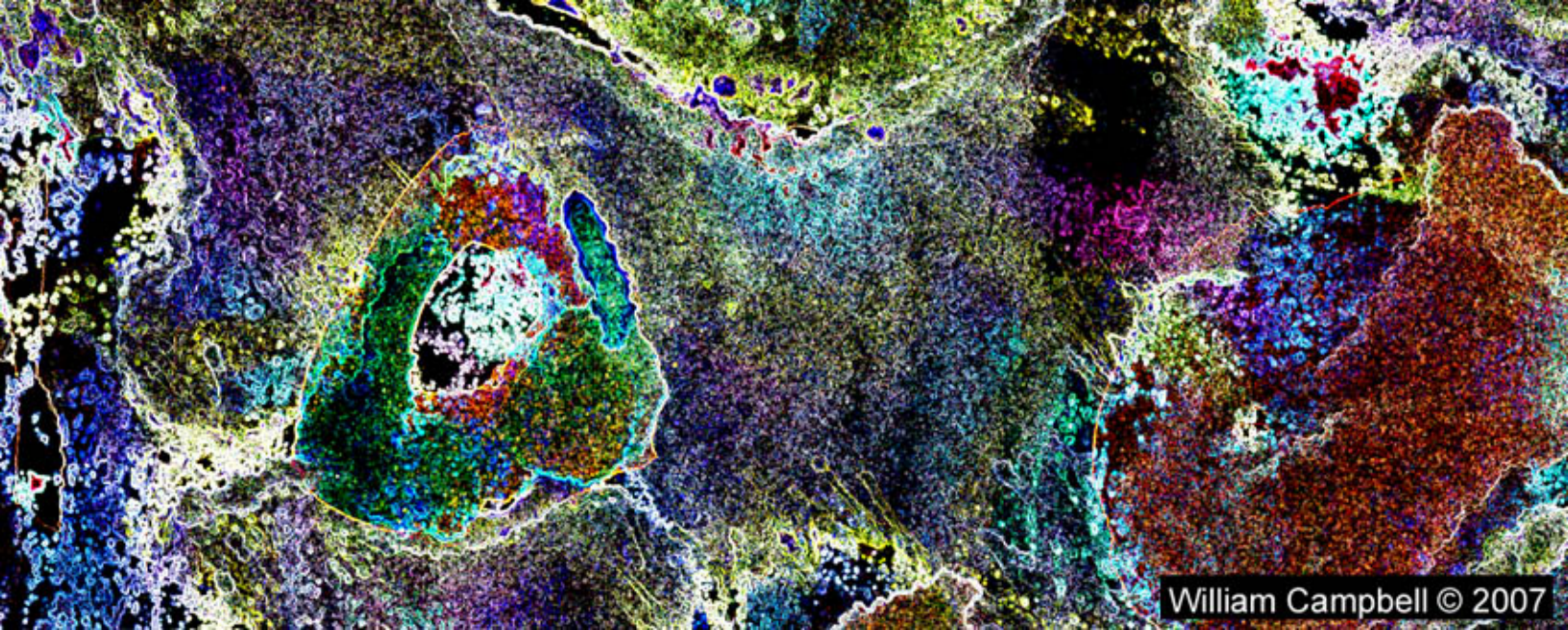




William Campbell © 2007