# User protocol for graphene FET GRFETS10

## **Product Info graphene FET S10**

Monolayer CVD grown Graphene based field effect transistors (FET) S10. Gate Oxide thickness: 90 nm Gate Oxide material: SiO<sub>2</sub> Resistivity of substrate: 1-10  $\Omega$ ·cm Metallization: Nickel/Aluminium 140 nm Graphene field-effect mobility: >1000 cm<sup>2</sup>/V·s Residual charge carrier density: <2 x 10<sup>12</sup> cm<sup>-2</sup> Dirac point: 10-40 V Yield >75 %

#### **Absolute Maximum Ratings**

Maximum gate-source voltage:  $\pm$  50 V Maximum temperature rating: 150 °C Maximum drain-source current density: 10<sup>7</sup> A/cm<sup>2</sup>

#### **Device configuration**

This Graphene FET chip provides 36 graphene devices distributed in a grid pattern on the chip. 30 devices have Hall-bar geometry and 6 have 2-probe geometry.

The Hall-bar devices can be used for Hall measurements as well as 4-probe and 2-probe measurements. There are graphene channels with varied dimensions to allow systematic investigation of device properties.





# **Channel geometries**



#### **Device Features**

- State-of-art graphene FETs utilizing consistent high-quality CVD grown monolayer graphene
- Devices are not encapsulated and can be functionalized by additives
- Perfect platform for sensor research and development
- 36 individual graphene FETs per chip
- Mobilities typically >  $1000 \text{ cm}^2/\text{V}\cdot\text{s}$

#### **Applications**

- Graphene device research<sup>1</sup>
- FET based sensor research for active materials deposited on graphene<sup>2</sup>
- Chemical sensors<sup>2</sup>
- Biosensors<sup>3,4</sup>
- Bioelectronics<sup>3,4</sup>
- Magnetic sensors
- Photodetectors

#### **Basic handling instructions**

The monolayer CVD graphene used in this FET device is highly prone to damage by external factors. To maintain the quality of the devices, we recommend taking the following precautions:

- Be careful when handling the graphene FET chip. Tweezers should not contact the device area directly.
- Metallic tweezers should be avoided, since they can damage/scratch the chip's surface
- Take precautions against electrostatic discharge

• Ideally store in inert atmosphere or under vacuum in order to minimize adsorption of unknown

- species from the ambient air
- Do not sonicate the graphene FET dies
- Do not apply any plasma treatment to the graphene FET dies
- Do not subject the graphene FET dies to strongly oxidizing reagents





## **Usage protocol:**

## 2-probe devices

These devices allow field-effect measurements by simultaneously applying two voltages:

 $\bullet$  Source-drain voltage (V\_{SD}): applied between the two probes (source and drain), while one of them is grounded (see Figure below).



Scheme of the 2-probe device, with the corresponding electrical measurement configuration.

 $V_{SD}$  enables the transport of charge carriers through the graphene channel, generating an associated source-drain current ( $I_{SD}$ ).  $V_{SD}$  can be varied in order to get the desired  $I_{SD}$  outcome (see Figure below).



Typical output curve measured at room temperature in vacuum.

• Gate voltage ( $V_G$ ): applied to the Si on the substrate.  $V_G$  creates an electric field on the graphene channel, modulating the conductivity of graphene (see Figure below).



Typical transfer curve measured at Vsp= 20mV (right), measured at room temperature in vacuum.

The Si can be contacted either from the top surface by scratching the 90 nm-thick  $SiO_2$  with a diamond pen in one of the chip corners; or alternatively from underneath the chip, for instance via a probe station chuck.

#### Hall Bars

#### 1. Field-effect measurement

A common improvement of the 2-probe GFET measurement is to apply a source-drain voltage between two outer contacts, measure the current between those two contacts, while measuring the voltage directly across the

graphene channel using two additional inner contacts,  $V_{12}$  (see Figure below). This way, the resistance of the graphene channel can be measured without inducing any voltage drops at the graphene-metal interfaces.





Scheme of the 4-probe measurement in a Hall bar device, with the corresponding electrical measurement configuration

The graphene-metal interface resistance does depend on  $V_G$  but not in the same way as the graphene channel resistance. Therefore measuring the graphene channel resistance directly in the 4-probe measurement configuration can achieve greater sensitivity to applied gate fields or surface charge changes.

The resistivity of graphene is usually expressed per unit thickness, i.e. the so-called sheet resistance:  $R_s = R_{CH} \frac{W}{L1}$ 

wherein  $R_{CH}$  is the resistance of the graphene channel, W and  $L_1$  is the width and inner length of the graphene channel, respectively. The field-effect mobility ( $\mu_{FE}$ ) can be calculated using the following equation:

$$\mu_{FE} = g \cdot \frac{1}{C_{SiO2}},$$

wherein:

- $g = d_{\sigma}/dV_{G}$  is the transconductance, with  $\sigma = 1/R_{S}$ ,
- $C_{SiO2}$  is the capacitance per unit area of the 90 nm-thick SiO<sub>2</sub> dielectric.

 $\mu_{12}$  is usually calculated using the maximum transconductance.

The field-effect charge carrier density (n\_{\rm FE}) is calculated as follows:  $n_{FE} = \mu_{FE} \cdot R_S/e$ 

In order to extract the residual carrier concentration  $n_0$ , i.e. the charge carrier density at the Dirac point, we can use the following equation:  $n_{FE} = \sqrt{n_0^2 + n_G^2}$ ,

wherein  $n_G$  is the gate-induced charge carrier density, which is calculated from: wherein  $V_D$  is the Dirac voltage and  $v_F$  is the Fermi velocity.

# 2. Hall Effect measurement

Hall Effect measurement is an alternative way to obtain the mobility and charge carrier density of graphene.

In this case,  $V_{SD}$  is applied between the horizontal contacts, whereas the transversal voltage or Hall voltage,  $V_H$ , is measured.  $V_H$  varies with the applied out-of-plane magnetic field B, due to the Lorentz force the charge carriers experienced. The Lorentz force deflects the charge carries toward the transverse contact, an electric field is thus created and measured by  $V_H$  (see below figure).



 $V_G - V_D = \frac{e}{C_{SiO2}} n_G + \frac{\hbar v_F \sqrt{\pi \cdot n_G}}{e},$ 



Scheme of the Hall Effect measurement configuration in a Hall bar.

 $n_H =$ The Hall mobility ( $\mu_H$ ) and charge carrier density ( $n_H$ ) are calculated according to:

$$R_H = \frac{dV_H}{dB} \cdot \frac{1}{I_{SD}}$$

Where  $R_H$  is the Hall coefficient:

$$\mu_H = \frac{n_H \cdot e}{R_s}$$

The mobility can then be calculated as:

# **De-Doping treatment**

Graphene on SiO<sub>2</sub> is often p-doped after exposing to air due to the adsorption of water molecules and other adsorbents that shifted the Dirac point to positive gate voltages, which can cause the Dirac voltage fall out of the recommended gate voltage range. Simultaneously, a large hysteresis can usually be observed between the forward and backward sweeps of the transfer curve.

Immersing the graphene FET chip in acetone for at least 12h can reduce doping. After that, the chip should be rinsed with IPA, dried with an Ar or  $N_2$  gun, prior to the measurement of the chip. In order to keep this treatment effective, electrical characterization should be carried out in inert atmosphere or in vacuum.

In addition, storage of the chips in a low humidity environment ( $N_2$  cabinet, desiccator, or vacuum) is highly recommended.

## **Reference:**

1. Wafer-scale statistical analysis of graphene field-effect transistors-part II: analysis of device properties By Smith, Anderson D.; Wagner, Stefan; Kataria, Satender; Malm, B. Gunnar; Lemme, Max C.; Ostling, Mikael From IEEE Transactions on Electron Devices (2017), 64(9), 3927-3933.

2. MoS2-graphene heterostructures as efficient organic compounds sensing 2D materials By Pham, Tung; Ramnani, Pankaj; Villarreal, Claudia C.; Lopez, Jhoann; Das, Protik; Lee, Ilkeun; Neupane, Mahesh R.; Rheem, Youngwoo; Mulchandani, Ashok From Carbon (2019), 142, 504-512.

3. All-electrical graphene DNA sensor array By Abbott, Jeffrey; Ham, Donhee; Xu, Guangyu From Methods in Molecular Biology (New York, NY, United States) (2017), 1572(Biosensors and Biodetection, Volume 2), 169-187.

4. Review-Field-Effect Transistor Biosensing: Devices and Clinical Applications By Syu, Yu-Cheng; Hsu, Wei-En; Lin, Chih-Ting From ECS Journal of Solid State Science and Technology (2018), 7(7), Q3196-Q3207.

